



# High Sensitivity Protein 250 Kit for 2100 Bioanalyzer Systems

## Kit Guide

*For Research Use Only - Not for use in diagnostic procedures*



# Notices

## Document Information

Document No: SD-UF0000063  
Rev. D.00  
Edition: 10/2023

## Copyright

© Agilent Technologies, Inc. 2000-2023

No part of this manual may be reproduced in any form or by any means (including electronic storage and retrieval or translation into a foreign language) without prior agreement and written consent from Agilent Technologies, Inc. as governed by United States and international copyright laws.

## Warranty

The material contained in this document is provided "as is," and is subject to being changed, without notice, in future editions. Further, to the maximum extent permitted by applicable law, Agilent disclaims all warranties, either express or implied, with regard to this manual and any information contained herein, including but not limited to the implied warranties of merchantability and fitness for a particular purpose. Agilent shall not be liable for errors or for incidental or consequential damages in connection with the furnishing, use, or performance of this document or of any information contained herein. Should Agilent and the user have a separate written agreement with warranty terms covering the material in this document that conflict with these terms, the warranty terms in the separate agreement shall control.

## Technology Licenses

The hardware and/or software described in this document are furnished under a license and may be used or copied only in accordance with the terms of such license.

## Restricted Rights Legend

U.S. Government Restricted Rights. Software and technical data rights granted to the federal government include only those rights customarily provided to end user customers. Agilent provides this customary commercial license in Software and technical data pursuant to FAR 12.211 (Technical Data) and 12.212 (Computer Software) and, for the Department of Defense, DFARS 252.227-7015 (Technical Data - Commercial Items) and DFARS 227.7202-3 (Rights in Commercial Computer Software or Computer Software Documentation).

## Safety Notices

### CAUTION

A **CAUTION** notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in damage to the product or loss of important data. Do not proceed beyond a **CAUTION** notice until the indicated conditions are fully understood and met.

### WARNING

A **WARNING** notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in personal injury or death. Do not proceed beyond a **WARNING** notice until the indicated conditions are fully understood and met.

# Contents

<b>1</b>	<b>High Sensitivity Protein 250 Kit</b>	<b>5</b>
<b>2</b>	<b>Required Equipment for High Sensitivity Protein 250 Assay</b>	<b>7</b>
<b>3</b>	<b>High Sensitivity Protein 250 Assay Principles</b>	<b>8</b>
<b>4</b>	<b>General workflow for the High Sensitivity Assay</b>	<b>9</b>
<b>5</b>	<b>Setting up Assay Equipment and Bioanalyzer</b>	<b>10</b>
	Setting up the Chip Priming Station	11
	Setting up the Bioanalyzer	12
	Starting the 2100 Expert Software	13
<b>6</b>	<b>Essential Measurement Practices</b>	<b>14</b>
<b>7</b>	<b>High Sensitivity Labeling Protocol</b>	<b>16</b>
	Preparation of Optimal Conditions for Labeling	16
	Verification of Absence of Interfering Substances	17
	Verification of Sample pH	18
	pH Adjustment for Samples	19
	Buffer Systems Suitable for the Labeling Reaction	22
	Labeling Reaction with Fluorescent Dye	23
<b>8</b>	<b>Protocol for On-Chip Analysis of Labeled Proteins</b>	<b>26</b>
	Dilution of Labeled Proteins	27
	Alternative Workflows	28
	Preparing the Denaturing Solution	29
	Preparing the Samples and the Ladder	30
	Loading the Gel Matrix	31
	Loading the Samples	33
	Inserting a Chip in the 2100 Bioanalyzer Instrument	34

Starting the Chip Run 35  
Cleaning Electrodes after a Chip Run 37

**9 Checking your High Sensitivity Protein 250 Results 38**

High Sensitivity Protein 250 Ladder Well Results 38  
High Sensitivity Protein 250 Sample Well Results 40  
Quantitative Evaluation of Sample Results 42

**10 Compatibility List for the Labeling Reaction 45**

**11 List of Known Effects from Dilution Buffers 47**

## 1

# High Sensitivity Protein 250 Kit

**Table 1 High Sensitivity Protein Kit**

<b>Agilent High Sensitivity Protein 250 Kit (5067-1575)</b>	
<b>High Sensitivity Protein Chips</b>	<b>Agilent High Sensitivity Protein 250 Labeling Reagents (5067-1577)</b>
10 Chips	● (green) 10x Protein 250 Standard Labeling Buffer (10xSLB), 10-fold concentrate (1 vial)
1 Electrode Cleaner	○ (clear) Ethanolamine (1 vial)
	● (blue) DMSO (1 vial)
<b>Syringe Kit</b>	● (blue) Labeling Dye (1 vial, separate light-tight bag)
1 Syringe	<b>High Sensitivity Protein 250 Ladder (5067-1578)</b>
	● (yellow) Ladder (1 vial, sufficient for three labeling reactions)
	<b>High Sensitivity Protein 250 Reagents (for On-Chip analysis, 5067-1576) &amp; Supplies</b>
	● (red) Gel-Matrix (1 vial, prefiltered)
	● (purple) Destaining Solution (1 vial)
	○ (white) Sample Buffer (3 vials)
<b>Prerequisites for labeling procedure</b>	
Type	Lysates, Extracts, Column Fractions, purified Proteins, lyophilized Proteins
Concentration	1 ng/μL to 3 μg/μL total protein
Volume	5 μL per labeling reaction
pH Value	Adjustment to pH 8.0 – 9.0
Optimal Matrix	Standard labeling buffer, supplied
Interferences	All components with primary amino or thiol groups

**Table 2 Physical Specifications**

Type	Specification
Analysis run time	30 min
Samples per chip	10
Sample volume	5 $\mu$ L
Kit stability	6 months
Kit size	10 chips 10 sample/chip = 100 samples/kit
Compatible buffers	"Compatibility List for the Labeling Reaction" on page 45

**Table 3 Specifications for analysis of labeled proteins**

Type	Specification
Sizing range	10 – 250 kDa
Typical sizing resolution	10 %
Typical sizing accuracy	10 % (BSA <sup>1</sup> )
Sizing precision	3 % CV (BSA)
Sensitivity (Signal/Noise > 3)	1 pg/ $\mu$ L (labeled BSA) in water on chip, 5 pg/ $\mu$ L (labeled BSA) in PBS on chip, labeling reaction at 1 ng/ $\mu$ L of total protein
Quantitative range	up to 4 orders of magnitude (0.1 – 1000 ng/ $\mu$ L BSA)
Quantitative precision	20 % CV (BSA)

<sup>1</sup> BSA = Bovine Serum Albumin

## 2

# Required Equipment for High Sensitivity Protein 250 Assay

Equipment Supplied with the Agilent 2100 Bioanalyzer System

- Chip priming station (5065-4401)

Additional Material Required (Not Supplied)

- 0.5 mL tubes
- Vortexer
- Deionized water
- Microcentrifuge
- 0.5 mL heating block or water bath (95 – 100 °C)
- pH-meter or indicator strips (basic range)

Further Information

Visit the Agilent website. It offers useful information, support, and current developments about the products and technology:

[www.agilent.com/en/product/automated-electrophoresis/bioanalyzer-systems](http://www.agilent.com/en/product/automated-electrophoresis/bioanalyzer-systems).

### 3

## High Sensitivity Protein 250 Assay Principles

The complete High Sensitivity Protein 250 kit contains chips and reagents for labeling of proteins with a dedicated fluorescent dye and subsequent sizing and quantitation.

The Assay workflow consists of two major steps:

- Covalent labeling of proteins with a fluorescent dye
- Separation and detection of labeled proteins with on-Chip-Electrophoresis

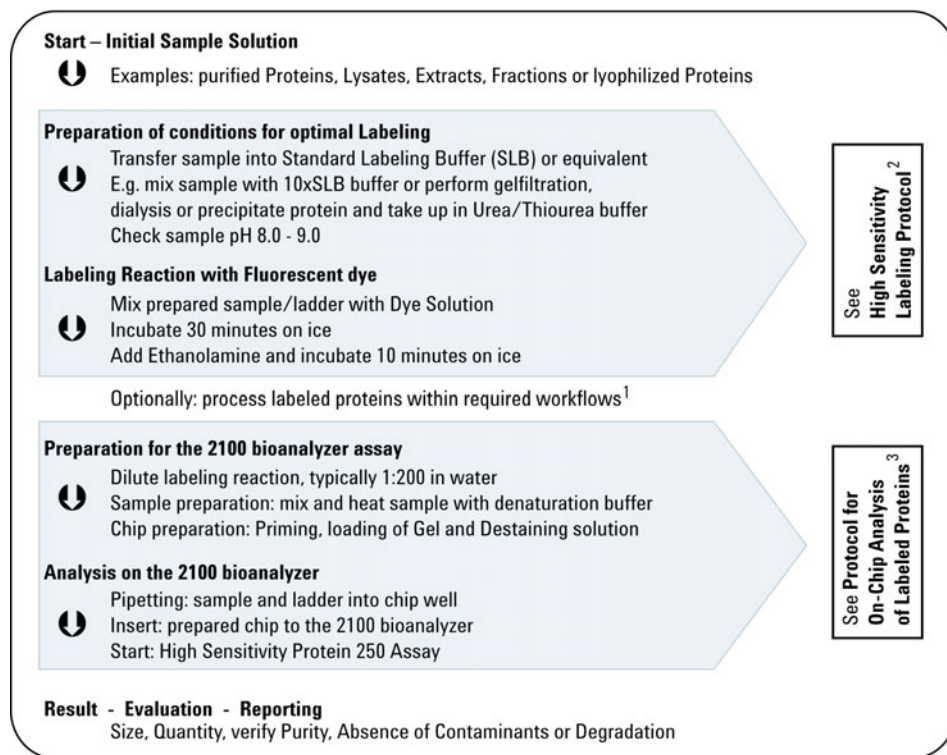
Epsilon-amino groups of lysine residues of proteins are covalently modified with a fluorescent dye by N-Hydroxy succinimidyl (NHS) ester chemistry.

Prepared 2100 Bioanalyzer chips consist of an interconnected set of polymer filled microchannels that sieve proteins by size as they are driven through it by means of electrophoresis. Agilent Protein kits are designed for use with the 2100 Bioanalyzer system only.



# 4

## General workflow for the High Sensitivity Assay



<sup>1</sup> Alternative workflows may be any purification, depletion or fractionation technique. See “Alternative Workflows” on page 28

<sup>2</sup> See “High Sensitivity Labeling Protocol” on page 16

<sup>3</sup> See “Protocol for On-Chip Analysis of Labeled Proteins” on page 26

## 5

# Setting up Assay Equipment and Bioanalyzer

Before beginning the labeling procedure:

You have to ensure that

- the sample is prepared for optimal labeling conditions (see “Preparation of Optimal Conditions for Labeling” on page 16).
- the labeling dye is reconstituted (see “Reconstitution of Fluorescent Dye with DMSO” on page 24).
- an ice bath is available

Before beginning the chip preparation protocol, ensure that the chip priming station and the Bioanalyzer are set up and are ready to use.

You have to

- replace the syringe at the chip priming station with each new protein kit
- adjust the base plate of the chip priming station to position A
- adjust the syringe clip at the chip priming station to the middle position
- adjust the Bioanalyzer's chip selector to position 1
- ensure that a heating block (95 – 100 °C) is available for heat denaturation
- start the 2100 Expert software (revision B.02.06 or higher), connect successfully to a 2100 Bioanalyzer and load the High Sensitivity Protein 250 assay before you load a chip.

### NOTE

The High Sensitivity Protein 250 assay is an assay which requires staining of proteins with a dedicated labeling protocol. The labeling is done prior to the analysis on chip. Please read this guide carefully and follow all instructions to guarantee satisfactory results.

## Setting up the Chip Priming Station

### NOTE

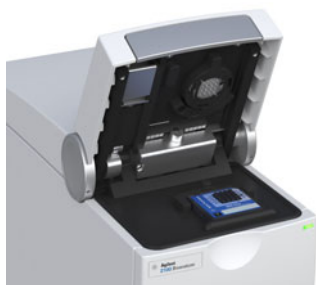
Replace the syringe with each new Reagent Kit.

- 1 Replace the syringe:
  - a Unscrew the old syringe from the lid of the Chip Priming Station.
  - b Release the old syringe from the clip. Discard the old syringe.
  - c Remove the plastic cap of the new syringe and insert it into the clip.
  - d Slide it into the hole of the luer lock adapter and screw it tightly to the Chip Priming Station.
- 2 Adjust the base-plate:
  - a Open the chip priming station by pulling the latch.
  - b Using a screwdriver, open the screw at the underside of the base plate.
  - c Lift the base plate and insert it again in position A. Retighten the screw.
- 3 Adjust the syringe clip:
  - a Release the lever of the clip and lift it up or down to adjust it to the middle position.



## Setting up the Bioanalyzer

- 1 Open the lid of the Bioanalyzer and make sure that the electrode cartridge is inserted in the instrument. If not, open the latch and insert the electrode cartridge.



**Figure 1** Electrode cartridge inserted in the instrument (graphic shows an example).

- 2 Remove any remaining chip.

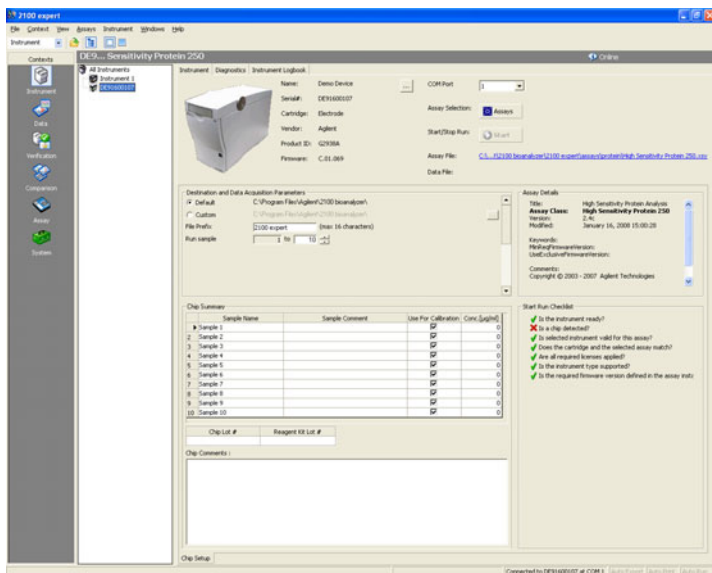
## Starting the 2100 Expert Software

To start the software:

- 1 Go to your desktop and double-click the following icon.



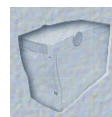
The screen of the software appears in the *Instrument* context. The icon in the upper part of the screen represents the current instrument/PC communication status:



Lid closed, no chip or chip empty



Lid open



Dimmed icon: no communication



Lid closed, chip inserted

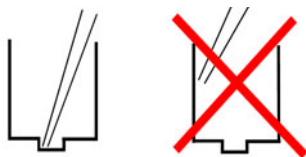
- 2 If more than one instrument is connected to your PC, select the instrument you want to use in the tree view.



## 6

# Essential Measurement Practices

- Handle and store all reagents according to the instructions under storage conditions.
- Avoid sources of dust or other contaminants. Foreign matter in reagents and samples or in the wells of the chip will interfere with assay results.
- Allow all reagents and samples to equilibrate to room temperature for 30 minutes before use and vortex.
- Always insert the pipette tip to the bottom of the chip well when dispensing the liquid. Placing the pipette at the edge of the well may lead to poor results.



- Use a new syringe and electrode cleaners with each new kit.
- Use loaded chips within 5 minutes. Reagents might evaporate, leading to poor results.
- Do not touch the 2100 Bioanalyzer instrument during analysis and never place it on a vibrating surface.
- Use 0.5 mL tubes to denature samples. Using larger tubes may lead to poor results, caused by evaporation.
- Keep suitable aliquots of the labeling reaction of the High Sensitivity Protein 250 Ladder undiluted at  $-28 - -15\text{ }^{\circ}\text{C}$  ( $-18 - 5\text{ }^{\circ}\text{F}$ ). Avoid freeze thaw cycles to prevent precipitation.
- The High Sensitivity Protein 250 Assay gel-matrix comes pre-filtered. It is ready to use after thawing.
- For Protein analysis under reducing conditions a 1 M DTT solution is required.
- Samples from labeling reactions need to be diluted prior to analysis. Do not dilute heat denatured samples.
- Relative concentrations indicated in the Peak Table of the Expert Software may need correction for the dilution step (e.g. 1:200). See [“Checking your High Sensitivity Protein 250 Results”](#) on page 38.
- Use calibrated pipets with proper pipetting technique. Choose suitable pipets and tips for low volumes.

Storage Conditions:

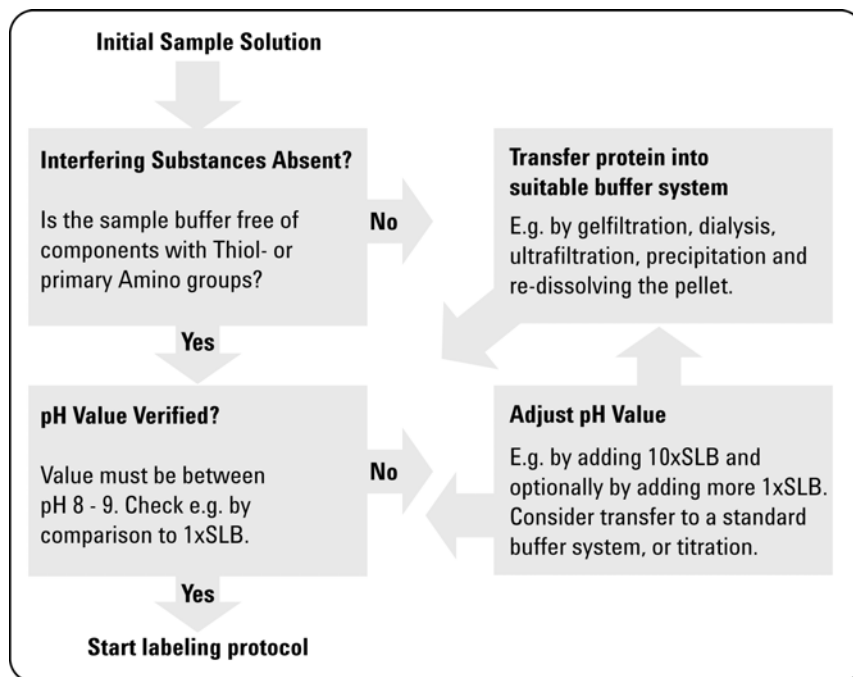
- Keep all reagents frozen at  $-28 - -15$  °C ( $-18 - 5$  °F) when not in use to avoid poor results caused by reagent decomposition.
- Avoid freeze thaw cycles for the ladder preparation.
- Protect all following reagents from light: labeling dye, sample buffer, destaining solution and any dye-labeled ladder/protein solution.
- Remove light covers only when pipetting. Dye decomposes when exposed to light.
- Store chips at room temperature.

# 7

## High Sensitivity Labeling Protocol

### Preparation of Optimal Conditions for Labeling

Follow all instructions for adjusting optimal labeling conditions when preparing the samples. It is crucial to have sample solutions in the pH range from pH 8 – 9 for satisfactory results of the complete assay. Note that it is essential as well to avoid interfering substances (see “Compatibility List for the Labeling Reaction” on page 45) in sample solutions.





## Verification of Absence of Interfering Substances

The solution with the sample protein should not contain disturbing components with Thiol- or primary Amino groups. Negative Interference on the labeling efficiency is known from primary amine or thiol groups of buffer components such as DTT,  $\beta$ -Mercaptoethanol, Glutathione, free amino acids and Imidazole. These buffer components compete with the sample protein in the reaction for the reactive labeling reagent (fluorescent dye). This competition lowers the effective reagent concentration and the amount of dye bound to protein, which subsequently lowers peak areas.

Interference from detergents such as CHAPS, Triton X-100, SDS or Tween 20 may have an influence the overall assay performance. Although labeling efficiency may be enhanced the analysis of labeled protein is affected in sizing and quantitation with on-Chip-Electrophoresis at higher detergent concentrations.

It is recommended to test any buffer deviating from the standard labeling buffer on its suitability for labeling by directly comparing it to a recommended buffer (see ["Buffer Systems Suitable for the Labeling Reaction"](#) on page 22).

### NOTE

It was found that a concentration of 1 mM DTT present in the Standard Labeling Buffer reduced the peak area of dedicated immunoglobulines by half. DTT contains thiol groups.

### NOTE

Find a comprehensive list of compatible sample buffers and possible effects on the labeling reaction under ["Compatibility List for the Labeling Reaction"](#) on page 45.

## Verification of Sample pH

Optimal labeling with the fluorescent dye will take place only if the pH is between pH 8.0 and pH 9.0. It is crucial to actively check the pH of sample solutions prior to the labeling reaction.

Verification of the pH value for the labeling reaction can be done by transferring a droplet to a pH-indicator strip which focuses on the basic range. Alternatively use a pH-meter on aliquots of your sample or buffer.

Labeling with the fluorescent dye will take place only if the pH is greater than pH 8.0. In pH-ranges beyond pH 9.0 other amino acid groups will be labeled as well. Both deviations from the optimum will negatively influence the assay results. Uniform labeling efficiency across all samples and the ladder is required for good quantitation results and reproducibility.

**NOTE**

For control of pH values special indicator strips focused on pH range 6.5 – 10.0 (0.2 – 0.5 unit resolution, Merck KGaA Darmstadt) were used successfully.

**NOTE**

Consider the usage of micro pH electrodes for measurements of small sample amounts.

**NOTE**

Labeling efficiency is greatly influenced by the pH value of the sample matrix.

## pH Adjustment for Samples

In case the sample shows a pH below pH 8.0 or above pH 9.0 it is necessary to adjust it to around pH 8.5 prior to the labeling reaction.

Dilution of sample with 10xSLB:

Standard labeling buffer is supplied as 10-fold concentrate (10xSLB, ●). To shift the pH add one part 10xSLB to 9 parts of sample. Consider this dilution factor in quantitation later on. Make sure the buffer capacity is suitable to shift the pH into a pH range between pH 8.0 and pH 9.0. Starting from the 10x concentrate the 1xSLB final concentration is 30 mM Tris. An estimation for the buffer capacities can be done from the molarities of the buffering substance in your individual sample buffer. The molarity of initial sample buffer substances should be much lower than 30 mM.

Confirm the adjusted pH prior to the labeling reaction by re-testing the solution. A good comparison for color pH-indicator strips is to test an aliquot of 1xSLB in parallel.

In case the pH adjustment with the above given recommendation is not applicable to your sample matrix consider to transfer sample proteins into a suitable buffer (see “[Buffer Systems Suitable for the Labeling Reaction](#)” on page 22).

**NOTE**

For example PBS (26mM  $\text{NaH}_2\text{PO}_4$ ; 41mM  $\text{Na}_2\text{HPO}_4$ , 79mM NaCl, pH 7.4) as sample matrix can be shifted in pH adequately by the addition of 1/10 volume 10xSLB to a pH > 8 which is suitable for labeling.

**NOTE**

Optionally more 1xSLB can be used to shift the pH furthermore. Do not exceed a final Tris concentration higher than 100 mM. Tris contains an amino group that may influence the labeling reaction at higher concentrations.

**NOTE**

In case addition of 10xSLB is not sufficient to shift the pH to the desired range, consider careful addition of 50 mM NaOH. Titrate your sample matrix offline to determine the required amount. Make sure the proteins under investigation are not harmed by this procedure.

**NOTE**

Deionized water should be at a neutral pH. Please check prior to use.

## Transfer of Protein into Suitable Buffer

In case the sample is in a buffer matrix that contains interfering substances it is recommended to transfer the proteins of interest to a suitable buffer, prior to the labeling reaction. Such suitable buffer may be the 1-fold concentrated standard labeling buffer (1xSLB), the recommended Sodiumbicarbonate buffer or the recommended Urea/Thiourea buffer (see “[Buffer Systems Suitable for the Labeling Reaction](#)” on page 22). Any other deviating buffer should be tested in direct comparison for equivalent efficiency in labeling.

Generally the following methods for transfer can be used:

### Gelfiltration

Gelfiltration e.g. in spin cartridges can be used for convenient removal of small sized interfering substances and buffer exchange. Equilibrate the resin with 1xSLB, Sodiumcarbonate or Urea/Thiourea buffer and proceed according to the manufacturers instructions. If buffer is exchanged completely to a recommended buffer, further pH adjustment is not necessary. Please verify the sample pH. Consider the final sample volume and dilution factors in quantitation.

### NOTE

The Zeba™ Micro Desalt Spin Columns were successfully used to transfer unknown samples into Standard Labeling Buffer or Sodiumbicarbonate buffer for the labeling reaction.

### Dialysis

Dialysis should be done with equilibrated membranes featuring suitable size cut-offs (size-retention e.g. at 3 kDa or 10 kDa) in accordance with the expected protein size. Dialysis should be performed for a sufficient time and with adequate buffer exchanges. If buffer is exchanged completely to a recommended buffer, further pH adjustment is not necessary. Please verify the sample pH.

### Ultrafiltration

Ultrafiltration e.g. in spin cartridges can be used for convenient re- concentration after a sample was diluted e.g. with 10x or 1x standard labeling buffer. Small interfering substance can be reduced in concentration by this procedure. The membrane should be equilibrated with e.g. 1xSLB and the cut-off (size-retention e.g. at 3 kDa or 10 kDa) should be in accordance with protein size.

## High Sensitivity Labeling Protocol

### Verification of Sample pH

Dissolving a pellet of sample proteins

This method requires precipitation and solubilization of sample proteins. Impurities are depleted effectively unless they are co-precipitated. Solubilization into the recommended Urea/Thiourea buffer (see “[Buffer Systems Suitable for the Labeling Reaction](#)” on page 22) can be done in small scale and enables concentration of the initial sample content. A simple acetone precipitation protocol is given below.

Simple clean-up protocol for proteins by acetone precipitation:

- Add 4-fold volume of chilled Acetone (-28 – -15 °C (-18 – 5 °F), not supplied)
- Incubate for ≥ 30 minutes at -28 – -15 °C (-18 – 5 °F)
- Pellet the protein by centrifugation: 10 minutes, 15000 g, 2 – 8 °C (36 – 46 °F)
- Optionally: Wash pellet with a chilled 4:1-mixture of Acetone:Water
- Remove supernatant completely and let pellet dry at room temperature
- Add Urea/Thiourea buffer to the pellet and dissolve to e.g. 1 µg/µL

Dissolving a lyophilisate

Take up a solid protein lyophilisate free of interfering substance into a suitable buffer e.g. 1xSLB. Lyophilization prerequisite is that the sample protein is dissolved in a completely volatile buffer. Verify absence of interfering substance in the lyophilisate. Note that not all proteins are stable during the freeze-drying process.

#### NOTE

The 2-D Clean-Up kit was successfully used to prepare unknown samples for the labeling reaction.

## Buffer Systems Suitable for the Labeling Reaction

Standard labeling buffer (SLB): 30 mM Tris/HCl, pH 8.5

The Standard labeling buffer is supplied with the kit as 10-fold concentrate (10xSLB; 300 mM Tris/HCl, pH > 8.5). It can be added to samples in order to shift the pH into the suitable range. To obtain a 1xSLB e.g. for buffer exchange in gel filtration or for comparison in pH-measurements, dilute the 10xSLB 1:10 with deionized water. E.g. add 10  $\mu$ L 10xSLB to 90  $\mu$ L water.

Sodiumbicarbonate buffer: 100 mM NaHCO<sub>3</sub>

Preparation of 50 mL: accurately weigh in solid Sodiumbicarbonate (NaHCO<sub>3</sub>, 420 mg), add water, stir until complete dissolution and fill up to 50 mL. Sodiumbicarbonate buffer decomposes with time. Prepare fresh frequently, store e.g. 2 weeks at 2 – 8 °C (36 – 46 °F) at the most. Check pH value to be between pH 8.0 and 9.0 prior to use.

Urea/Thiourea buffer: 30 mM Tris/HCl, 7 M Urea, 2 M Thiourea, pH 8.5

Preparation of 500  $\mu$ L: accurately weigh in solid Urea (210 mg) and Thiourea (76 mg), add 50  $\mu$ L of 10xSLB and add 250  $\mu$ L deionized water. Mix until complete dissolution of solid components and fill up with deionized water to 500  $\mu$ L. Urea/Thiourea buffer decomposes with time. Prepare fresh frequently, store e.g. 2 month at -28 – -15 °C (-18 – 5 °F) at the most. Check pH value to be between pH 8.0 and 9.0 prior to use.

---

**NOTE**

Deionized water should be at a neutral pH. Please check prior to use.

---

**NOTE**

Ammonium carbonate buffers are not suitable for the labeling reaction since the Ammonium competes with the protein for the reactive Fluorescent Dye and thus interferes with the labeling reactions.

---

## Labeling Reaction with Fluorescent Dye

### WARNING

#### Handling Reagents

The dye can cause serious eye damage. Because the dye binds to nucleic acids, it should be treated as a potential mutagen.

The ladder can cause eye irritation.

Kit components contain DMSO. DMSO is skin-permeable and can elevate the permeability of other substances through the skin.

Thiourea is harmful if swallowed, suspected of causing cancer and of damaging the unborn child, and is toxic to aquatic life with long lasting effects.

- ✓ Follow the appropriate safety procedures and wear personal protective equipment including protective gloves and clothes as well as eye protection.
  - ✓ Follow good laboratory practices when preparing and handling reagents and samples.
  - ✓ Always use reagents with appropriate care.
  - ✓ For more information, refer to the material safety data sheet (MSDS) on [www.agilent.com](http://www.agilent.com).
- 

## Labeling Reaction with Fluorescent Dye

Epsilon-amino groups of lysine residues of proteins are covalently modified with a fluorescent dye by N-Hydroxy succinimidyl (NHS) ester chemistry.

The Fluorescent dye reagent needs reconstitution in DMSO. For the labeling reaction it is mixed with sample solution or ladder. Excess dye will be quenched after the reaction by Ethanolamine. Ladder volume provided in this kit is sufficient for 3 independent labeling reactions. Each Ladder labeling reaction may be aliquoted and used successively for several chips.

## Reconstitution of Fluorescent Dye with DMSO

- 1 Remove the fluorescent dye and DMSO from the freezer (●, both vials).
- 2 Equilibrate both vials to room temperature, DMSO should be completely thawed.
- 3 Centrifuge dye vial at 10000 g for 2 min to collect solid dye particles at the bottom of the vial.
- 4 Visually localize the dye pellet and add 54  $\mu$ L DMSO onto the pellet.
- 5 Vigorously vortex the Dye solution until all solid components are completely dissolved.
- 6 Label the vial with date.
- 7 Store Dye solution in the dark at  $-28 - -15$  °C ( $-18 - 5$  °F) for up to 6 months, thaw completely prior to use.

### NOTE

DMSO is solid at  $2 - 8$  °C ( $36 - 46$  °F). Please thaw DMSO and dye solution completely. Vortex the reconstituted dye prior to use.

## Performing the Labeling Reaction

- 1 Thaw reconstituted dye solution completely and vortex prior to use.
- 2 Prepare 5  $\mu$ L of Ladder (●) in a tube.
- 3 Prepare 5  $\mu$ L of protein sample (ready for labeling, see “Preparation of Optimal Conditions for Labeling” on page 16) per tube.
- 4 Place tubes on ice.
- 5 Start the labeling reaction by adding 0.5  $\mu$ L of reconstituted dye solution (●) to tubes with samples or ladder, vortex and spin down for 5 seconds.
- 6 Incubate 30 minutes on ice.
- 7 In order to quench any excess of dye add 0.5  $\mu$ L of Ethanolamine (○), vortex and spin down for 5 seconds.
- 8 Incubate 10 minutes on ice.  
Labeling of ladder and protein sample is finalized.
- 9 Start analysis of the labeled ladder and store remaining solution aliquoted (1  $\mu$ L) at  $-28 - -15$  °C ( $-18 - 5$  °F).
- 10 Start analysis of the labeled products or store labeling reaction mixture at  $-28 - -15$  °C ( $-18 - 5$  °F).



## High Sensitivity Labeling Protocol

### Labeling Reaction with Fluorescent Dye

**NOTE**

Performing the reaction on ice ensures uniform and reproducible protein labeling.

---

**NOTE**

Consider the use of tubes with low binding capacity for work with minimal amounts of protein. Eppendorf Safe-Lock Protein LoBind 0.5 mL Microcentrifuge tubes were tested successfully for the labeling reaction, dilution steps and storage.

---

## 8

# Protocol for On-Chip Analysis of Labeled Proteins

The on-chip analysis of proteins labeled with Fluorescent Dye requires dilution of the initial labeling reaction and heat denaturation prior to sample loading to the prepared chip.

- 1 After completing the initial steps in “[Setting up Assay Equipment and Bioanalyzer](#)” on page 10, you can perform the steps, as described in the following procedures.
- 2 Allow the Protein 250 gel matrix and destaining solution to equilibrate to room temperature for 30 minutes. Both solutions need no further preparation prior to use.
- 3 Allow the High Sensitivity Protein 250 denaturing solution to equilibrate to room temperature for 30 minutes. Prior to the first usage a preparation is needed, as described below.

### NOTE

All kit components can be stored at -28 – -15 °C (-18 – 5 °F).

### NOTE

It is important that all the reagents have room temperature before starting the next steps.

### NOTE

Protect the dye containing solutions such as labeled samples, sample buffer, denaturing solution and destaining solution from light.

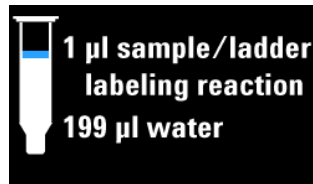
### NOTE

An entry of sample names and comments (e.g. dilution factor) is possible in the instrument context prior to starting a run.

## Dilution of Labeled Proteins

For analysis of samples directly from the labeling reaction:

- 1 Allow labeling reaction mix from the samples to equilibrate to room temperature.
- 2 Allow an aliquot from the ladder labeling reaction to equilibrate to room temperature.
- 3 Dilute sample and ladder 1:200 in water, e.g. add 1  $\mu\text{L}$  of the labeling reaction to 199  $\mu\text{L}$  water, vortex, and continue immediately with the on-chip analysis.

**NOTE**

In case a different dilution factor is considered in your customized workflow, please regard potential consequences (see [“High Sensitivity Protein 250 Sample Well Results”](#) on page 40). For alternative dilution buffers see [“List of Known Effects from Dilution Buffers”](#) on page 47.

**NOTE**

Dilution is necessary to avoid signal saturation and subsequent bias. Often, this is due to the Lower Marker peak, representing co-migrating excess dye from the labeling reaction and Lower Marker from the sample buffer.

**NOTE**

Diluted labeled sample and ladder should be analyzed immediately. Do not use this preparation after storage since proteins may precipitate, degrade or aggregate.

## Alternative Workflows

In case of proceeding alternative workflows with the labeled sample proteins the dilution step can be adapted. An alternative workflow may be any preparative technique. Labeling of lysates, extracts or other mixtures of proteins can also be done prior to their separation. Initially labeled proteins can serve as starting material in size exclusion, hydrophobic interaction, ion exchange chromatography, precipitation or separation according to isoelectric points. The multiple fractions of these approaches can subsequently be analyzed directly with the High Sensitivity Protein 250 Assay.

The described workflows usually implicate a depletion of the non-reacted dye. Therefore no or only little dilution is necessary.

## Preparing the Denaturing Solution

- 1 Remove one of the original vials of the High Sensitivity Protein 250 sample buffer (white ◯) or a vial with an aliquot of the High Sensitivity Protein 250 sample buffer from the freezer. Allow to equilibrate to room temperature for 30 minutes, then vortex.
- 2 For *reducing* conditions: To the 100  $\mu$ L sample buffer in the original vial (white ◯) add 3.5  $\mu$ L of 1 M Dithiothreitol (DTT) solution or 3.5 Vol.-% of 1 M DTT to an aliquot of sample buffer.

OR

For *non-reducing* conditions: Add deionized water to sample buffer instead of DTT.

- 3 Vortex for 5 seconds.



### NOTE

The denaturing solution contains a Fluorescent Dye functioning as lower marker in the assay. Protect this solution from light.

### NOTE

100  $\mu$ L denaturing solution is sufficient for 50 sample preparations. Store prepared denaturing solution at  $-28 - -15$  °C ( $-18 - 5$  °F).

### NOTE

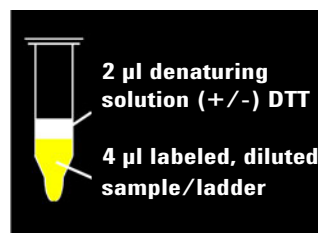
Avoid exposure of DTT containing buffer to the air since oxidation may degrade the reducing reagent leading to insufficient reduction of your sample proteins.

## Preparing the Samples and the Ladder

**Prerequisites**

For a list of compatible buffers for the dilution step or any matrix remaining from alternative workflows, please refer to “List of Known Effects from Dilution Buffers” on page 47.

- 1 Allow the denaturing solution (prepared as described above) to equilibrate to room temperature for 30 min, and vortex before use.
- 2 Combine 4  $\mu\text{L}$  of your labeled, diluted protein sample and 2  $\mu\text{L}$  of denaturing solution in a 0.5 mL microcentrifuge tube. Prepare multiples of this for replicates.
- 3 Combine 4  $\mu\text{L}$  of your labeled, diluted High Sensitivity Protein 250 Ladder and 2  $\mu\text{L}$  of denaturing solution in a 0.5 mL microcentrifuge tube.
- 4 Mix well and spin down for 15 seconds.
- 5 Place each sample tube and the ladder tube for 5 minutes in a heating block at 95 – 100 °C or in boiling water.



Ensure that the tubes are properly placed and heated. Do not heat for more than 5 minutes otherwise excessive evaporation might occur. The samples and ladder should not dry down.

- 6 Let the tubes cool down to ambient temperature.
- 7 Spin vials for 15 seconds to recover condensate or liquid.  
Samples and Ladder are now ready for chip loading.

**NOTE**

You will need 6  $\mu\text{L}$  sample preparation per chip well. No further dilution step is done. Prepare duplicates of samples or ladder in order to fill all 10 wells of a complete chip.

**NOTE**

Ladder performance is uniform for reducing and non-reducing conditions.

**NOTE**

Incubation time and temperature for heat denaturation may require optimization for your protein of interest.

## Loading the Gel Matrix

**NOTE**

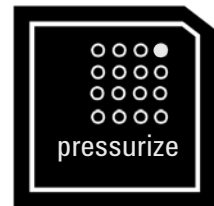
Before loading the gel matrix, make sure that the base plate of the chip priming station is in position (A) and the adjustable clip is set to the middle position. Refer to “Setting up the Chip Priming Station” on page 11 for details.

- 1 Allow the gel matrix to equilibrate to room temperature for 30 minutes before use. The gel comes prefiltered and is ready to use after equilibration to room temperature.
- 2 Take a High Sensitivity Protein chip out of its sealed bag and place it on the Chip Priming Station.
- 3 Pipette 12  $\mu$ L of the gel at the bottom of the well marked (●).

**NOTE**

When pipetting the gel matrix insert the tip of the pipette to the bottom of the chip well when dispensing. This prevents a large air bubble forming under the gel. Placing the pipette at the edge of the well may lead to poor results.

- 4 Set the timer to 90 seconds, make sure that the plunger is positioned at 1 mL and then close the Chip Priming Station. The lock of the latch will click when the priming station is closed correctly.
- 5 Press the plunger of the syringe down until it is held by the clip.
- 6 Wait for exactly 90 seconds and then release the plunger with the clip release mechanism.


**NOTE**

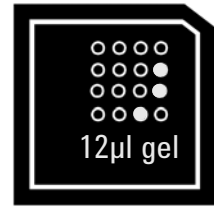
Use a new syringe with each new chip box. Mount syringe to the chip priming station and test this with the “seal test”. See details in the 2100 Expert Help Menu.


- 7 Visually inspect that the plunger moves back at least to the 0.3 mL mark.
- 8 Wait for 5 s, then slowly pull back the plunger to the 1 mL position.
- 9 Open the Chip Priming Station.

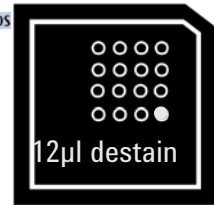
## Protocol for On-Chip Analysis of Labeled Proteins

### Loading the Gel Matrix

**10** Pipette 12  $\mu\text{L}$  of the gel mix in each of the 3 wells marked with .



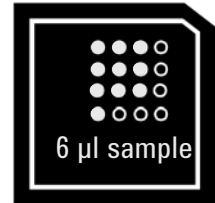
**11** Pipette 12  $\mu\text{L}$  of the destaining solution in the well marked .





## Loading the Samples

- 1 Pipette the complete volume of the heat denatured samples (6  $\mu$ L, prepared as described in “Preparing the Samples and the Ladder” on page 30) into the sample wells marked 1...10.




### NOTE

Complete volume from heat denaturation is potentially less than 6  $\mu$ L due to evaporation or pipetting inaccuracy. Use tight vials and calibrated pipets.

### NOTE

In case insufficient liquid levels are filled to the chip wells the 2100 Expert software will not start the chip run but give a warning. A loss of 5 % volume is acceptable.

- 2 Pipette the complete volume (6  $\mu$ L) of the heat denatured ladder (as described in “Preparing the Samples and the Ladder” on page 30) into the well marked with the ladder symbol .
- 3 Make sure that the run is started within 5 minutes. Select the High Sensitivity Protein 250 assay within the instrument context. Refer to the next topic on how to insert the chip in the 2100 Bioanalyzer instrument.



### NOTE

Do not leave any wells empty or the chip will not run. Pipette a sample or ladder replicate in any empty sample well.

## Inserting a Chip in the 2100 Bioanalyzer Instrument

- 1 Open the lid of the 2100 Bioanalyzer instrument.
- 2 Check that the electrode cartridge is inserted properly. Refer to “Setting up the Bioanalyzer” on page 12 for details.
- 3 Place the chip carefully into the receptacle. The chip fits only one way.

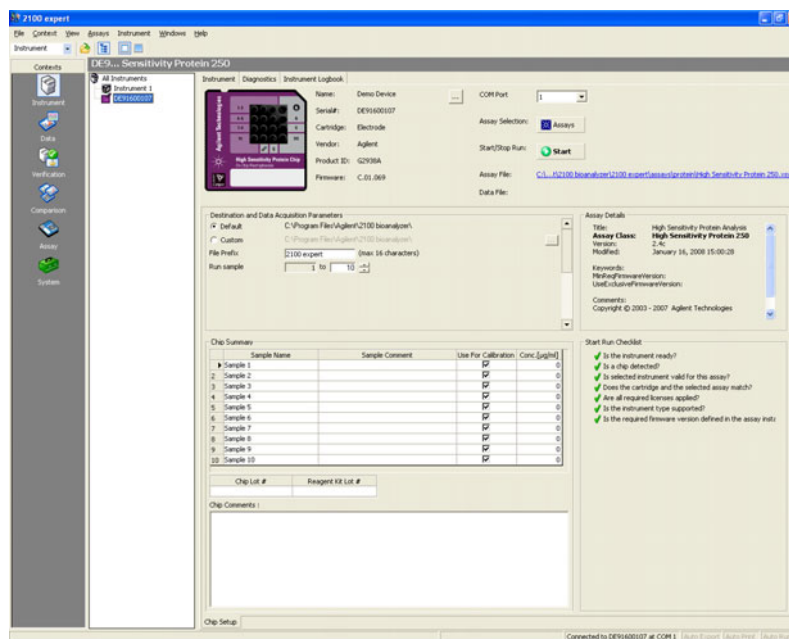
### CAUTION

#### Sensitive electrodes and liquid spills

Forced closing of the lid may damage the electrodes and dropping the lid may cause liquid spills resulting in bad results.

- ✓ Do not use force to close the lid and do not drop the lid onto the inserted chip.

- 4 Carefully close the lid. The electrodes in the cartridge fit into the wells of the chip.
- 5 The 2100 Expert software screen shows that you have inserted a chip and closed the lid by displaying the chip icon at the top left of the *Instrument* context.

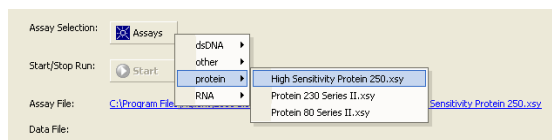


## Starting the Chip Run

### NOTE

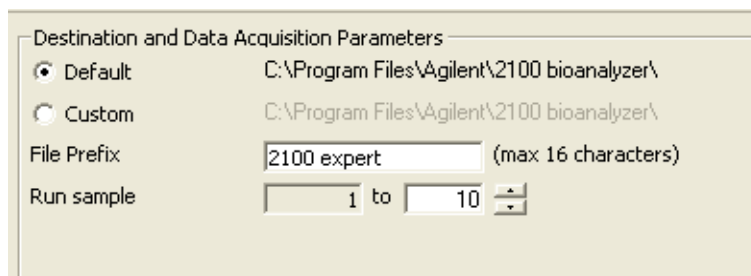
Please note that the order of executing the chip run may change if the Agilent Security Pack software (only applicable for 2100 Expert software revision B.02.06 and higher) is installed. For more details please read the 'User's Guide' which is part of the Online Help of your 2100 Expert software.

- 1 In the **Instrument** context, select the High Sensitivity Protein 250 assay from the Assay menu.



- 2 Accept the current **File Prefix** or modify it.

Data will be saved automatically to a file with a name using the prefix you have just entered. At this time you can also customize the file storage location and the number of samples that will be analyzed.



- 3 Click the **Start** button in the upper right of the window to start the chip run. The incoming raw signals are displayed in the **Instrument** context.



- 4 To enter sample information like sample names and comments, select the *Data File* link that is highlighted in blue or go to the **Data** context and select the **Chip Summary** tab. Complete the sample name table.

### NOTE

An entry of sample names and comments is also possible in the instrument context prior to starting a run. The information (e.g. dilution factor as comment) can be entered in the chip summary tab while the analysis takes place as well or after the run is finished.

**NOTE**

For aspects of absolute quantitation please refer to “Absolute quantitation” on page 43. For information on adjustment of the ladder concentration refer to “Adjustment of the ladder concentration” on page 44.

- 5 To review the raw signal trace, return to the *Instrument* context.

**CAUTION****Contamination of electrodes**

Leaving the chip for a period longer than 1 hour (e.g. over night) in the Bioanalyzer may cause contamination of the electrodes.

- ✓ Immediately remove the chip after a run.

- 6 After the chip run is finished, remove the chip from the receptacle of the Bioanalyzer and dispose of it according to good laboratory practices.

## Cleaning Electrodes after a Chip Run

When the assay is complete, *immediately* remove the used chip from the 2100 Bioanalyzer instrument and dispose of it according to good laboratory practice. Then perform the following procedure to ensure that the electrodes are clean (no residues are left over from the previous assay).

### NOTE

Use a new electrode cleaner with each new kit.

### CAUTION

#### Leak currents between electrodes

Liquid spill may cause leak currents between the electrodes.

✓ **Never fill more than 350  $\mu$ L of water in the electrode cleaner.**

- 1 Slowly fill one of the wells of the electrode cleaner with 350  $\mu$ L deionized analysis-grade water.
- 2 Open the lid and place the electrode cleaner in the 2100 Bioanalyzer instrument.
- 3 Close the lid and leave it closed for about 10 seconds.
- 4 Open the lid and remove the electrode cleaner.
- 5 Wait another 10 seconds to allow the water on the electrodes to evaporate before closing the lid.

### NOTE

After each chip run, empty and refill the electrode cleaner.

After 10 chip runs, replace the used electrode cleaner by a new one.

### NOTE

When switching between different assays, a more thorough cleaning may be required. Refer to the maintenance chapter and Troubleshooting Guide for details. It is part of the Online Help of the 2100 Bioanalyzer software.

# 9

## Checking your High Sensitivity Protein 250 Results

### High Sensitivity Protein 250 Ladder Well Results

To check the results of your run, select the Gel or Electropherogram tab in the Data context. The electropherogram of the ladder well window should resemble the one shown below.

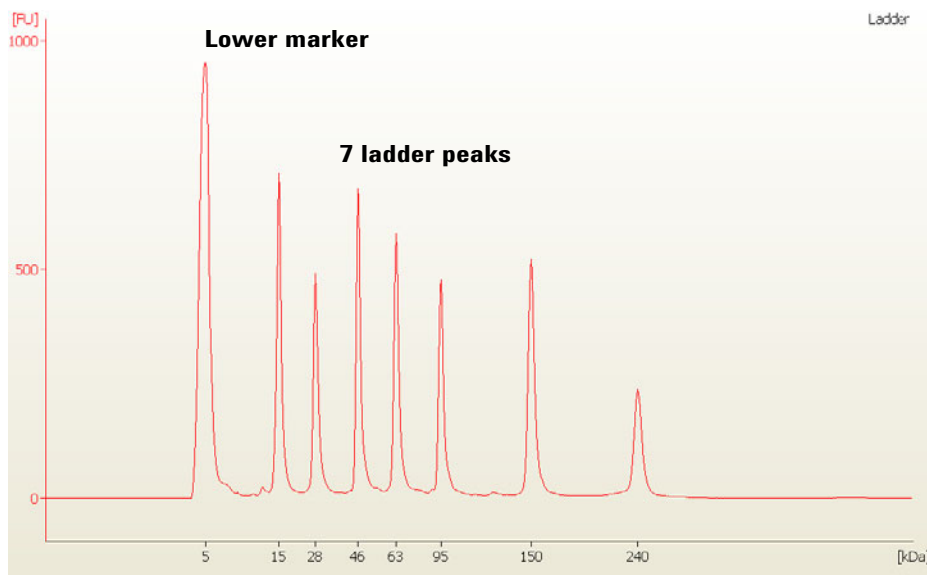


Figure 2 High Sensitivity Protein 250 ladder

## Checking your High Sensitivity Protein 250 Results

### High Sensitivity Protein 250 Ladder Well Results

Major features of a successful ladder run are:

- 7 ladder peaks and all peaks are well resolved
- 1 lower marker peak, usually higher than the ladder peaks
- Flat baseline
- Readings at 50 units higher than baseline readings

If the electropherogram of the ladder run does not resemble the one shown above check the handling of the labeling reaction and the handling of the on-chip analysis. For general assistance refer to the *2100 Expert Maintenance and Troubleshooting Guide* within the online Help of the 2100 Expert software.

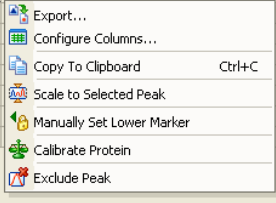
#### NOTE

The High Sensitivity Protein 250 assay uses the ladder for relative quantification. It works without upper marker as internal standard. Alignment of Ladder and Samples is based on the Lower marker only.

In case wrong peaks are identified as ladder peaks, exclude them by doing the following:

- 1 Move the cursor over the peak in the peak table and click the right mouse button.
- 2 Select *Exclude Peak* from ladder to make the change come into effect.

	Size [kDa]	Observations	
1	5.0	Lower Marker	
	9.4		
3	15.0	Ladder Peak	
4	28.0	Ladder Peak	
5	46.0	Ladder Peak	
6	63.0	Ladder Peak	
7	95.0	Ladder Peak	
8	150.0	Ladder Peak	
9	240.0	Ladder Peak	



## High Sensitivity Protein 250 Sample Well Results

To review the results of a specific sample, select the sample name in the tree view and highlight the Results sub-tab. The electropherogram of the sample run should resemble the one shown below for the High Sensitivity Protein 250 assay.

Major features of a successful protein sample run are:

- A lower marker peak is visible.
- Baseline prior to the lower marker peak is flat.
- All sample peaks are migrating later than the lower marker peak and are resolved from it (depending on sample).
- The lower marker migrates between 16 and 24s (analysis turned off).
- The baseline reading is between 5 and 80 FU (analysis turned off).

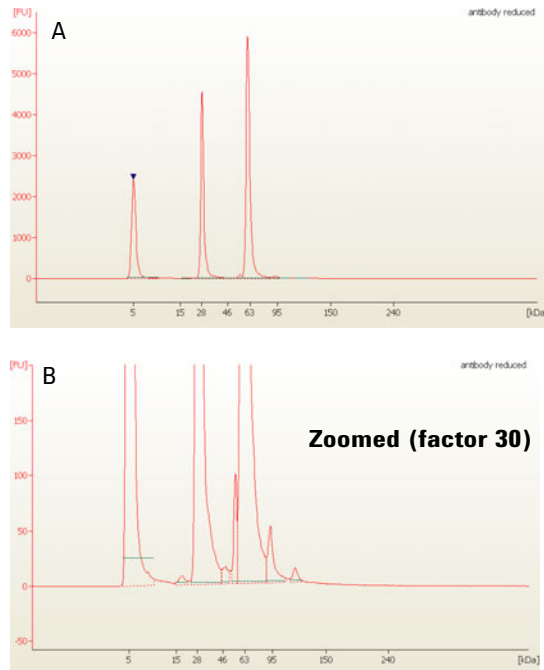


Figure 3 Sample lane, reduced antibody (A) full scale (B) Zoom-In

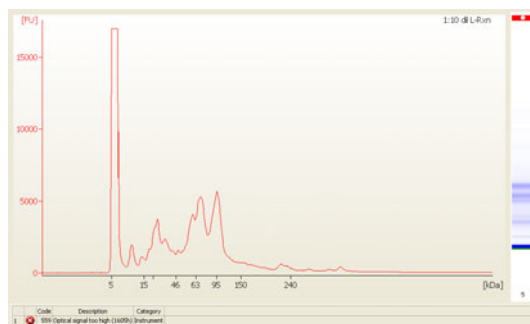


**NOTE**

The Lower marker peak represents lower marker (a fluorescent dye) in the sample buffer plus excess of labeling dye. Therefore the overall lower marker peak height is higher in case of low protein input to the labeling reaction compared to reactions with high protein input.

Insufficient dilution of labeling reactions will yield a saturation of the 2100 Bioanalyzer fluorescence detector and subsequently an error message (*optical signal too high*). The electropherogram will show rectangular peak shapes for the lower marker or for dominant sample peaks. Sizing of all peaks will be affected since peak finding for the lower marker is biased. Quantitation of distorted sample peaks is affected due to the peak shape. Tailing of the lower marker peak beyond 10 kDa may disturb integration of smaller protein peaks.

Dilution of labeling reaction by 1:200 is recommended (page 24). For alternative workflows it must be estimated or determined experimentally.



**Figure 4** Detector saturation, choose higher dilution factor

**NOTE**

In case the Lower marker peak disturbs the analysis excess of labeling dye from the reaction mixture may be depleted. Gel filtration can be employed to yield smaller lower marker peaks and to avoid the saturation of the fluorescence detector. This approach generally allows application of lower dilution factors and may increase sensitivity.

## Quantitative Evaluation of Sample Results

### Relative quantitation

For the High Sensitivity Protein 250 assay quantitation is done with the help of the ladder time corrected area. The sum for the identified 7 ladder peaks is compared with the sample peak. The area under the “lower” marker is not taken into consideration. Because the concentration of the ladder is known (4167 pg/μL; initial ladder concentration is  $10^6$  pg/μL and undergoes 5/6 dilution during labeling and 1:200 in dilution with water), the concentration for each sample peak can be calculated by the rule of three.

The result table gives the relative concentration. By default it refers to the concentration of the sample solution after dilutions. Please multiply the result with the dilution factors (standard protocol: 5/6 *labeling reaction* plus 1:200 *water dilution* equals 1:240 in total) to receive the concentration of your initial sample solution.

Besides a relative quantitation based on the ladder, an absolute quantitation is available using external standard proteins (see “[Absolute quantitation](#)” on page 43).

Internal standards spiked into the sample solution prior to the labeling reaction enable an alternative quantitation strategy. Introduction of a known peak concentration and employing the “% of Total” result from the analysis allows concentration calculation. The approach using an internal standard allows to control for matrix-related labeling effects.

#### NOTE

Correct assignment of the ladder peaks is required for correct relative quantitation. Please check this for each run.

## Checking your High Sensitivity Protein 250 Results

### High Sensitivity Protein 250 Sample Well Results

#### Absolute quantitation

If absolute quantitation is required with a standard protein, mark the check box **Use For Calibration** and enter standard concentration.

	Sample Name	Sample Comment	Use For Calibration	Conc. [µg/ml]	Status	Observation	Result Label	Result Color
1	IgG 1000 pg/ul	standard; dilution 1:200	<input checked="" type="checkbox"/>	1000	✓		Contains heavy + light chain	
2	IgG 100 pg/ul	standard; dilution 1:200	<input checked="" type="checkbox"/>	100	✓		Contains heavy + light chain	
3	IgG 10 pg/ul	standard; dilution 1:200	<input checked="" type="checkbox"/>	10	✓		Contains heavy + light chain	
4	IgG 1 pg/ul	standard; dilution 1:200	<input checked="" type="checkbox"/>	1	✓		Contains heavy + light chain	
▶	blank	control; dilution 1:200	<input type="checkbox"/>	0	✓		no peaks at 25+50 kDa	
6	sample IgG non-reduced	unknown; dilution 1:200	<input type="checkbox"/>	0	✓		Contains heavy + light chain	
7	sample IgG reduced	unknown; dilution 1:200	<input type="checkbox"/>	0	✓		Contains heavy + light chain	
8	sample IgG reduced	unknown; dilution 1:200	<input type="checkbox"/>	0	✓		Contains heavy + light chain	
9	sample IgG non-reduced	unknown; dilution 1:200	<input type="checkbox"/>	0	✓		Contains heavy + light chain	
10	blank	control; dilution 1:200	<input type="checkbox"/>	0	✓		no peaks at 25+50 kDa	

Chip Lot #  Reagent Kit Lot #

Chip Comments :

Sample Information | Study Information | Instrument Information | Standard Curve | Calibration Curve

Import... Export...

Local | **Global**

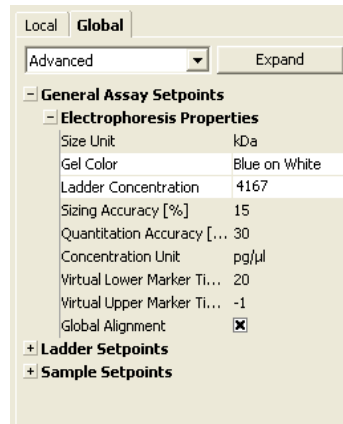
Advanced  Expand

- + General Assay Setpoints
- + Ladder Setpoints
- Sample Setpoints
  - + Alignment
  - + Quantitation
  - + Sizing
  - Calibration
    - Calibration - Expected ... 0
    - Calibration - Peak Time -1
    - Calibrate All
  - + Smear Analysis
  - + Baseline Calculation
  - + Filter Settings
  - + Baseline Correction
  - + Integrator

The highest peak in the calibration samples will be used for a linear calibration curve, which will be available on a new tab in the Chip Summary. The linear regression from time corrected areas of the calibration peaks are applied to the time corrected area of any sample peak to calculate the peak concentration.

Peaks from samples showing the same size as the calibration peak or manually assigned peaks will be quantified using the calibration curve. By default only one peak per lane is recognized as calibrated protein. In case all sample peaks should be evaluated with the calibration curve select the function "calibrate all" under advanced global setpoints from the setpoint explorer.

### Adjustment of the ladder concentration



The dilution step of labeling mixes of sample and ladder is variable by the nature of the assay. The proposed dilution is 1:200 in water. Preset ladder concentration in the 2100 Expert software is thus 4167 pg/μL. If different dilution ratios are used the total ladder peak concentration may be adjusted accordingly under advanced setpoints.

Another helpful customization is to adjust the ladder concentration setpoint to  $10^6$  pg/μL, i.e. the initial ladder concentration prior to the labeling reaction. Manual correction for the dilution factors (standard protocol: *5/6 labeling reaction* plus *1:200 water dilution* equals 1:240 in total) thus can be omitted. The concentrations given by the 2100 Expert Software will now refer to the initial sample concentration prior to the labeling reaction as long as samples are treated exactly like the ladder.

The Ladder Concentration setpoint may be considered as an adjustable factor for calibration purpose as well. In comparison to a well characterized master sample the ladder concentration value can be adjusted to let the assay yield the assumed concentration of this master sample. This adjusted value can be re-used in subsequent runs if the identical labeled ladder preparation is applied in the ladder well. This procedure allows to recycle a calibration result over several individual chips. Consider to generate a larger amount of labeled ladder in one reaction, e.g. 15 μL, to yield a sufficient number of identical aliquots.

## Compatibility List for the Labeling Reaction

The following table lists protein sample buffers and buffer components which are known to have low impact on the labeling reaction. Others have a medium or strong negative influence.

Please refer to the preparation of recommended buffers, see “Buffer Systems Suitable for the Labeling Reaction” on page 22.

**Table 4** Compatibility List for the Labeling Reaction

Low Impact: 50% - 150% signal compared to 1xSLB <sup>1</sup>
30 mM Tris-HCl, pH 8.5 (Standard Labeling Buffer, 1xSLB <sup>*</sup> ).
100 mM Sodiumbicarbonate <sup>*</sup> pH 9.0
1xSLB, 7 M Urea, 2 M Thiourea (Urea/Thiourea Buffer <sup>*</sup> )
1xSLB with 0.1 mM DTT
1xSLB with 1.0 mM DTT
1xSLB with 20 mM EDTA
1xSLB, 0.04% Sodiumazide
1xSLB, 1% CHAPS
1xSLB, 1% Triton
1xSLB, 1% Tween
100 mM Tris-HCl, pH 8.5
1xSLB + PBS (30 mM Tris-HCl, 26 mM NaH <sub>2</sub> PO <sub>4</sub> , 41 mM Na <sub>2</sub> HPO <sub>4</sub> , 79 mM NaCl) pH 8.5
50 mM HEPES, pH 8.0
50 mM CHES, pH 8.5
30 mM Tris-HCl, 1.25 M NaCl, pH 9.0
30 mM Tris-HCl, 30% Glycerol, pH 8.3
30 mM Tris-HCl, 0.9 M KCl, pH 9
30 mM Tris-HCl, 50 mM MgCl <sub>2</sub> , pH 8.3

## Compatibility List for the Labeling Reaction

### High Sensitivity Protein 250 Sample Well Results

**Table 4** Compatibility List for the Labeling Reaction

**Low Impact: 50% - 150% signal compared to 1xSLB<sup>1</sup>**

*Medium Impact: 10% - 50% signal compared to 1xSLB\**

1xSLB with 10 mM DTT

50 mM Tris-HCl, 100 mM NaCl, 20 mM Glutathion, pH 8.5

1xSLB, 1% SDS, pH 8.7

30 mM Tris-HCl, pH 7.3

*Strong Impact: < 10% signal compared to 1xSLB\**

100 mM Glycine/NaOH, pH 9

50 mM Tris-HCl, 500 mM NaCl, 500 mM Imidazole, pH 9

<sup>1</sup> These are recommended buffers for the labeling reaction.

## List of Known Effects from Dilution Buffers

The following table lists tested matrices for on-chip analysis of labeled proteins, i.e. the compatible buffers for dilution of the labeling reaction or the matrix after following an alternative workflow.

**Table 5** Tested matrices

Dilution Buffer	Sizing Effect	Comment
Water <sup>1</sup> , de-ionized	0%	Reference
<b>Low Impact, Signal Intensity &gt; 50%</b>		
10 mM Tris, 1 mM EDTA (TE)	< 3%	
10% DMSO, water	< 3%	
50 mM CHES, pH 8.5	< 3%	
4 M Urea, water	< 3%	
40% Acetonitrile, 0.1% formic acid	< 3%	
1% SDS in water	< 3%	
30% Glycerol in water	< 3%	
<b>Medium Impact, Signal Intensity 10% - 50%</b>		
1 mg/ml BSA <sup>2</sup> , PBS (26 mM NaH <sub>2</sub> PO <sub>4</sub> , 41 mM Na <sub>2</sub> HPO <sub>4</sub> , 79 mM NaCl) pH 7.4	< 3%	
1xSLB	< 3%	
1xSLB, 100 mM NaCl	< 3%	
1xSLB, 100 mM KCl	< 3%	
1xSLB, 100 mM DTT	< 3%	
100 mM Sodiumbicarbonate pH 9	< 3%	
PBS	< 3%	
PBS, 250 mM Imidazole	< 3%	
50 mM Na-Acetate, pH 5	< 10%	LM migration affected

## List of Known Effects from Dilution Buffers

### High Sensitivity Protein 250 Sample Well Results

**Table 5**    **Tested matrices**

Dilution Buffer	Sizing Effect	Comment
1xSLB, 20 mM EDTA	< 10%	
PBS, 0.1% Tween-20	< 10%	
50 mM HEPES pH 7.5	< 10%	LM migration affected
50 mM MES, pH 6	< 10%	LM migration affected
50 mM Sodium-Phosphate, pH 7.5	> 10%	LM migration affected
50 mM MOPS, pH 7	> 10%	Migration time and Peak shape affected
<b>Strong Impact, Signal Intensity &lt; 10%</b>		
50 mM Tris-HCl, 100 mM NaCl, 20 mM Glutathion, pH 8	< 3%	High ionic strength affects sample injection
30 mM Tris-HCl pH 8.5, 50 mM MgCl <sub>2</sub>	< 3%	High ionic strength affects sample injection

<sup>1</sup> This is the recommended solution for the dilution.

<sup>2</sup> Non-labeled proteins will not be detected in the assay.



## In This Book

you find the procedures to analyze Protein samples with the Agilent High Sensitivity Protein 250 Kit and the Agilent 2100 bioanalyzer.

- Agilent High Sensitivity Protein 250 Kit
- Assay Principles
- General workflow
- Labeling Protocol
- On-Chip Analysis of Labeled Proteins
- Agilent High Sensitivity Protein 250 Results
- Compatibility List for the Labeling Reaction
- List of Known Effects from Dilution Buffers

[www.agilent.com](http://www.agilent.com)

© Agilent Technologies Inc. 2000-2023  
Edition: 10/2023

Document No: SD-UF0000063 Rev. D.00

