



NOTICE: ProZyme was purchased by Agilent in July 2018. Documents for products and product lots manufactured before August 2019 will contain references to ProZyme. For more information about these products and support, go to: www.agilent.com/en/contact-us.



Rapid N-Glycan Preparation with 2-AB

Enzymatic deglycosylation, fluorescent labeling with 2-AB (2-aminobenzamide) and cleanup of excess dye for analysis by LC and other methods.

- Optional sample purification for monoclonal antibodies and Fc-fusion proteins
- Non-selective, rapid release and recovery of N-Glycans from up to 96 glycoprotein samples at a time using a microplate centrifuge
- Optimized reaction conditions help to preserve labile glycans, such as sialic acid
- Non-selective chemistry for stoichiometric labeling of glycans, independent of structure
- Complete labeling in 1 hour using Rapid-Reductive-Amination™
- Purified 2-AB-labeled N-glycans are eluted in water, ready for analysis
- Flexible, high-throughput format: process 1 to 192 samples per run (2 Kits simultaneously)
- Compatible with microplate liquid handling on a broad range of automation platforms

Product Code: GP24NG-AB

TABLE OF CONTENTS

| | page |
|--|------|
| Kit Contents. | 2 |
| Storage Requirements | |
| Additional Required Reagents/Equipment | |
| Optional Reagents and Supplies | |
| Safety and Handling. | 4 |
| Introduction. | 5 |
| Using the Kit. | 6 |
| Preparation of Samples | |
| Enzyme Incubation | |
| Labeling Incubation | |
| Protocols. | 10 |
| Overview of the Procedure | |
| Getting Started | |
| Purify (optional) | |
| Digest | |
| Label | |
| Cleanup | |
| Analysis of Labeled N-Glycans. | 32 |
| Tips and Hints. | 32 |
| References. | 34 |
| Technical Assistance. | 35 |
| License to Use. | 36 |
| Trademarks and Tradenames. | 37 |
| Product Use and Warranty. | 38 |
| Other ProZyme Products & Kits. | 38 |
| Ordering Information. | 38 |

This product is intended for in vitro research use only.

NOTE: The following suggestions and data are based on information we believe to be reliable. They are offered in good faith, but without guarantee, as conditions and methods of use of our products are beyond our control. We recommend that the prospective user determine the suitability of our materials and suggestions before adopting them on a commercial scale. Suggestions for use of our products or the inclusion of descriptive material from patents and the citation of specific patents in this publication should not be understood as recommending the use of our products in violation of any patent or as permission to license to use any patents of ProZyme, Inc.

KIT CONTENTS

NOTE: We want successful results for our customers, so please read this entire booklet before starting the procedure.

| Item | Qty |
|---|----------|
| G5524-60001 Kit AssayMAP PA50 (purchased separately) | Optional |
| G5524-60001 AssayMAP PA50 (24 Cartridges) | |
| WS0296-24 Protein A Solution Set | |
| WS0251-24 Eluent (7.5 ml) | |
| WS0294-24 10x Wash Buffer (4 ml) | |
| GP24NG-AB GlykoPrep® Rapid N-Glycan Preparation with 2-AB | |
| GS24-RX GlykoPrep Digestion Module | 1 ea |
| WS0283 Digestion (RX) Cartridges (2 x 12 Cartridges) | |
| WS0256-24 Immobilization Reagent Set | |
| WS0226-24 Denaturation Reagent (8 ml) | |
| WS0255-24 Blocking Reagent (1.5 ml) | |
| WS0259-24 Digestion Reagent Set | |
| WS0278-24 N-Glycanase® (75 µl) | |
| WS0276 25x Digestion Buffer (700 µl) | |
| WS0229 Finishing Reagent (1 ml) | |
| Aluminum Sealing Film (1) | |
| GS24-AB GlykoPrep Rapid-Reductive-Amination 2-AB Labeling Module | 1 ea |
| WS0302-24 2-AB Solution (75 µl) | |
| WS0300-24 Reductant Solution (75 µl) | |
| GS24-CU GlykoPrep Cleanup Module | 1 ea |
| WS0293 Cleanup (CU) Cartridges (24 Cartridges) | |
| Aluminum Sealing Film (1) | |

Storage Requirements

This Kit is a mixed-temperature shipment (2–30°C). Upon arrival, store components as indicated. For best results, equilibrate materials to ambient temperature prior to use. The 2-AB Solution and Reductant Solution are hygroscopic and the 2-AB Solution is light-sensitive; please store these reagents at -20°C in their original packaging.

Additional Required Reagents/Equipment

Heater and Incubation Blocks, capable of 50–100°C, available from ProZyme as Product Code GS150
AssayMAP Labware: Racks, Receiver Plates and Lids
Other Labware: Waste Plates/Cleanup Collection Plates, 450 µl well volume (Thermo Fisher Scientific part number 07-202-502/Corning part number 3343 or equivalent)
Gilson Diamond® D200 Pipet Tips

NOTE: AssayMAP Labware to use with this Kit is available from ProZyme as a complete Starter Set (Product Code AM200), or purchased separately in sets of 10.

Centrifuge (capable of 50–1000 x g) and deep microplate rotor with a height clearance of ≥44 mm
Ultrapure, deionized water (Milli-Q® or equivalent)
Acetonitrile (100%, HPLC-grade)
Pipettors & disposable tips (P5/P10, P200 and P1000)
Vortexer or plate shaker
Nitrile gloves
Vial, 1-ml, polypropylene for use with organic solvents
Fume hood (for Labeling procedure)
Miscellaneous labware for buffers and dilutions
Glass, graduated cylinder, 25 ml

Microplate-compatible, centrifugal evaporator (e.g., SpeedVac® or similar) for drying finished N-glycans prior to labeling. A rotor that holds two shallow-well microplates and fits into a small-capacity rotary evaporator may be purchased from Thermo Fisher Scientific (Savant Model RH2MP).

Optional Reagents and Supplies

Multichannel pipettors & disposable tips (P5/P10 and P200) (Gilson or equivalent, compatible with Gilson Diamond® D200 pipette tips)
Microplate reader (capable of reading A₂₈₀) for measurement of antibody concentration after Purification
Pipette basins (must be polypropylene for use with organic solvents)

SAFETY AND HANDLING

Some of the reagents in this Kit are hazardous. Please refer to the Safety Data Sheets (SDS) included with the Kit or posted on ProZyme's website under the component name or Product Code.

<http://www.prozyme.com>

General Laboratory Procedures

Use powder-free, nitrile gloves for all sample handling procedures. Ensure that all glass, plasticware and solvents are free of glycosidases and environmental carbohydrates.

All procedures involving Labeling Reagents (2-AB Solution, Reductant Solution and 2-AB Labeling Reagent) should be performed in a dry environment with dry glassware and plasticware, using appropriate personal safety protection, eyeglasses and nitrile gloves, and where appropriate, in a fume hood.

INTRODUCTION

The GlykoPrep Sample Preparation Platform (GlykoPrep) dramatically streamlines glycoanalysis by facilitating optional protein purification, quantitative deglycosylation and separation of N-glycans, complete fluorescent labeling and efficient cleanup to reduce excess reagent peaks.

GlykoPrep is modular and can be integrated into any workflow, regardless of throughput or sample type. In order to match any standard sample preparation, Kit components are also available individually as the AssayMAP PA50 (for purification of Fc-containing antibodies only), Digestion Module and dye-specific Labeling & Cleanup Modules.

GlykoPrep is built on AssayMAP technology, microchromatography in a 96-well format, capable of automated high throughput. GlykoPrep may be performed using centrifugation to move liquid through the Cartridges (spin format), or with the Syringe Head on the Agilent AssayMAP Bravo Liquid Handling Workstation (GlykoPrep-plus). Using the spin format with a microplate centrifuge, up to 192 samples can be processed simultaneously with two 96-count Kits.

Important general information for achieving success with the spin format, as well as special tips particular to individual Modules, may be found in the GlykoPrep Guidebook under Using Specific Kits and Modules:

<http://www.prozyme.com/documents/TNGP100.pdf>

We also provide a modified Microfuge Method useful for those interested in using the spin format to run only a handful of samples with a benchtop microfuge and a PCR heater:

http://www.prozyme.com/documents/2-AB_Microfuge_Method.pdf

USING THE KIT

GlykoPrep Rapid N-Glycan Preparation with 2-AB combines the Digestion Module, the Rapid-Reduction-Amination 2-AB Labeling Module and the Cleanup Module, which may be purchased individually. The Labeling and Cleanup Modules may be purchased together as GP24-ABCU. Optional purification modules may be employed just prior to Digestion to allow glycoanalysis directly from cell culture as a single workflow (directions included for your convenience). For information on purification modules under development, please contact us.

Preparation of Samples

Sample Quantities

The quantitative binding for each Cartridge is:

| | |
|----------------------------|------------------------------------|
| AssayMAP PA50 Cartridge | 125 µg of MAb or Fc-fusion protein |
| RX Cartridge | 50 µg of most standard proteins |
| CU Cartridge | 30 µg of N-glycans |

NOTE: The binding capacity for specific glycoproteins may need to be determined.

Cartridges are capable of binding more target, but will do so with increasing breakthrough, making the process non-quantitative.

For quantitative loading, prepare an excess of 10% or more sample, and prepare replicates together. For example, for Digestion, samples should be denatured together and loaded individually.

Less than the maximum quantity may be processed, for example, when the sample is available only in limited amounts. The smallest amount of sample that will give good results depends on the sensitivity requirements of the analytical methods and the specific application (e.g., screening vs. QC release).

Sample Denaturation

Prior to deglycosylation, the samples are denatured by pre-mixing with Denaturation Reagent. The suggested sample concentration prior to deglycosylation is 1–5 mg/ml, and sufficient reagents have been provided for the standard sample concentration range.

NOTE: If quantitation is desired, pipetting less than 10 µl is not recommended; pipetting smaller volumes introduces variability, especially when samples are highly concentrated. If necessary, dilute the sample to within the 1-5 mg/ml range with Digestion Buffer before starting.

The Kit is useful for very dilute samples without requiring further concentration, by expanding this load step to multiple spins. See the GlykoPrep Guidebook section “Loading.”

When performed in a single spin, the amount loaded to each RX Cartridge should be 10–100 µl. The recommended starting ratio of Denaturation Reagent to sample is 1:1 (v/v).

NOTE: The final denatured Sample must be at least 50% Denaturation Reagent.

Example 1:
Sample concentration 1 mg/ml
Sample amount needed: 50 µg

50 µl (50 µg) Sample + 50 µl Denaturation Reagent
= 100 µl denatured sample

Example 2:

Sample concentration 5 mg/ml

Sample amount needed: 50 µg

10 µl (50 µg) Sample + 90 µl Denaturation Reagent

= 100 µl denatured sample

The current protocol employs a 5-minute, relatively gentle denaturation, but any custom denaturation may be performed and the subsequent protocol followed as described, as long as no SDS or other detergents are used. Please see the GlykoPrep Guidebook under Digestion Modules or contact us to discuss custom denaturation conditions for your glycoprotein.

Enzyme Incubation

Time

The Digest procedure has been optimized to deliver deglycosylation of N-glycans in 15-60 minutes. The optimal incubation time will vary depending on the specific glycoprotein; those which have proven to be resistant to deglycosylation via conventional enzymatic methods may require longer incubation times (up to 60 minutes). For glycoproteins that are comparatively easy to deglycosylate, such as monoclonal antibodies, a 15-minute incubation is generally sufficient. The selected Incubation Time will be used in the Digestion Module.

NOTE: It is critical not to exceed a 60-minute incubation, as the Cartridge resin bed may dry out, yielding uncertain results.

Finishing Reagent converts the glycosylamine produced by N-Glycanase digestion to a free reducing end, required for labeling with 2-AB via Rapid-Reductive-Amination. The

incubation time for Finishing is fixed at ten minutes.

Temperature

The GS150 Heater and Incubation Blocks are specially designed to provide rapid heat transfer through the Receiver Plate and into the packed bed of each Cartridge. The Incubation Blocks are sold separately (ProZyme Product Code WS0272) and can be used in any standard dry-block heater of the proper size, or pre-heated and used in an oven. Custom Incubation Blocks compatible with robotic systems are also available from ProZyme.

NOTE: If using the Microfuge Method format, PCR heat block substitutes for the GS150 Heater (with Incubation Blocks).

The GS150 Heater (with Incubation Blocks) is set to 50°C for the Digest procedure (deglycosylation and Finishing step). Please allow a minimum of 1 hour to equilibrate the Blocks before use. The Incubation Blocks have been designed with a thermometer well in the corner. We have verified that when the thermometer reads 50°C, the temperature in the Cartridge is ~37°C, the optimal temperature for both deglycosylation and in the well for finishing. If using a different heater, confirm the block temperature.

Labeling Incubation

Time

The Labeling incubation time is fixed at 60 minutes.

Temperature

The Labeling procedure requires the Incubation Blocks to be equilibrated at 70°C. If using a different heater, check the temperature. We have verified that when the block is at 70°C, the temperature in the well is ~54°C, the optimal temperature for 2-AB labeling.

PROTOCOLS

Overview of the Procedure

Purify (optional, purchased separately)

Antibodies or Fc-fusion proteins may be purified from crude samples using Protein A.

Digest

Samples (antibodies or other proteins) are denatured and immobilized.

N-glycans are released by N-Glycanase[®] digestion and eluted, finished and dried prior to labeling.

Label

Dried N-glycans are labeled with 2-AB by Rapid-Reductive-Amination.

Cleanup

Buffer salts and excess labeling reagents are removed; labeled N-glycans are eluted in water, ready for analysis.

Getting Started

Heater Setting

Turn on the GS150 Heater (with Incubation Block). Set to 50°C and allow to equilibrate for a minimum of 1 hour.

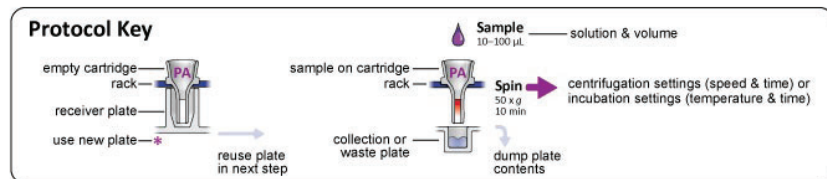
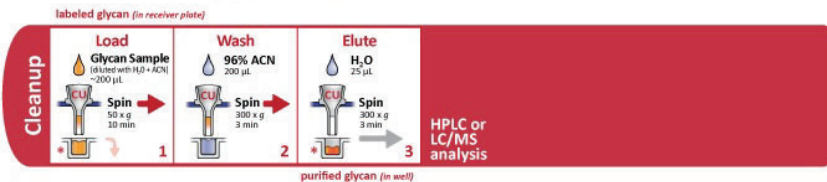
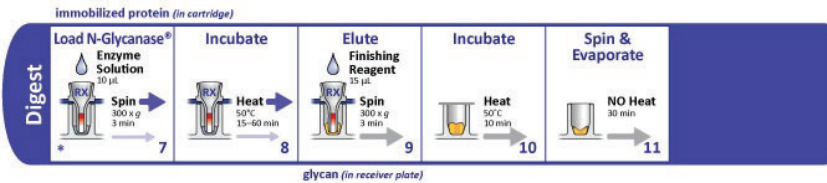
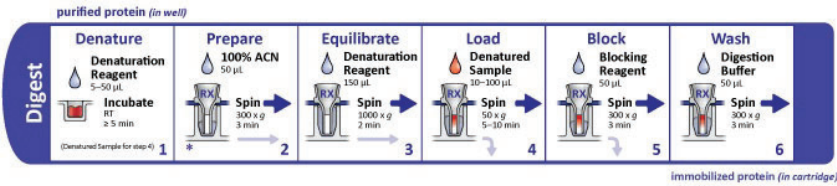
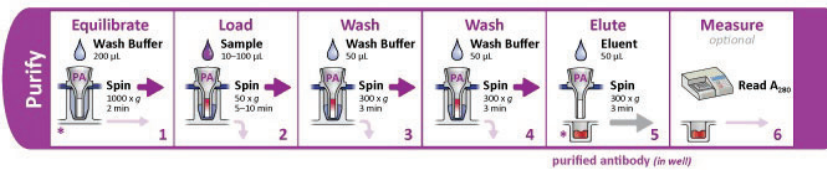
Centrifuge Settings

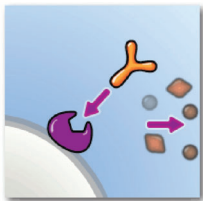
If the centrifuge does not have x g settings, determine the setting and the specific microplate rotor combination by consulting the operation manual or the manufacturer's website:

_____ rpm = 50 x g

_____ rpm = 300 x g

_____ rpm = 1000 x g



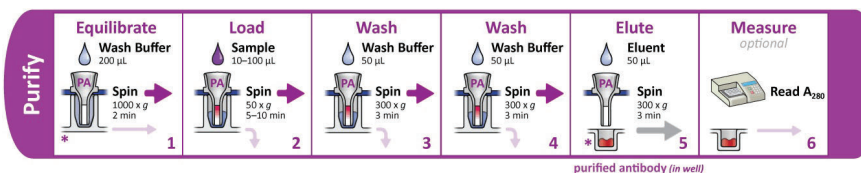


Purify (optional)

Protein A purifies antibodies or Fc-fusion proteins from cell-culture supernatants. All other samples must be purified by other methods; proceed to Digestion Module.

Overview

- 1 Equilibrate
- 2 Load
- 3 Wash
- 4 Wash (second time)
- 5 Elute
- 6 Measure (optional)



Reagents and other Supplies

AssayMAP PA50 Cartridges (supplied in the G5524-60010 Kit, 1 per sample)

Prepare two balanced AssayMAP PA50 Cartridge assemblies (Cartridges on Racks on Receiver Plates with Lids)

Wash Buffer (prepared below)

Eluent (supplied with the Kit)

NOTE: The prepared eluent must NOT contain glycine because it leaves a signature peak on the LC chromatogram.

Purification Collection Plate (supplied in the AM200 Starter Labware Set, or equivalent)

UV-compatible, flat-bottom, half-area plate for direct protein assay (optional)

PCR plate (optional)

Crude antibody or Fc-fusion protein samples (samples may contain NO MORE than 125 µg total protein, the quantitative binding capacity of the AssayMAP PA50 Cartridge; the amount loaded onto the downstream RX Cartridge may contain NO MORE than 50 µg total protein, the quantitative binding capacity of the downstream RX Cartridge). Samples should be between pH 6.5 and 8.5 and clear of particulates.

Preparation of Reagents

Wash Buffer

NOTE: May be prepared up to one week before use. Store at 2–8°C.

10x Wash Buffer (supplied with the Kit)
Ultrapure water

Dilute one volume of 10x Wash Buffer stock with nine volumes of ultrapure water to obtain Wash Buffer. Specifically, add 4 ml of 10x Wash Buffer stock to 36 ml of ultrapure water to make 40 ml of Wash Buffer.

For fewer samples, prepare 400 µl of Wash Buffer for each sample to be processed.

Procedure

Equilibrate

- 1.a Pipet 200 μl of Wash Buffer into the Sample Cup of each AssayMAP PA50 Cartridge.
- 1.b Spin at 1000 x g for 2 minutes; do not empty the Receiver Plates.

Load

- 2.a Load 10–100 μl of sample into the Sample Cup of each AssayMAP PA50 Cartridge (see Sample Loading Technique in the GlykoPrep Guidebook).
- 2.b Remove the Racks from the Receiver Plates. Empty the Receiver Plates and blot with a paper towel to avoid cross-contamination. Replace Racks.
- 2.c Spin at 50 x g until all Sample Cups are empty. The estimated spin time is 5 minutes for volumes between 10 and 50 μl or 10 minutes for volumes up to 100 μl .

Wash

- 3.a Pipet 50 μl of Wash Buffer into the Sample Cup of each AssayMAP PA50 Cartridge.
- 3.b Empty the Receiver Plate and blot with a paper towel.
- 3.c Spin at 300 x g for 3 minutes.

Wash (second time)

- 4.a Pipet 50 μl of Wash Buffer into the Sample Cup of each AssayMAP PA50 Cartridge.
- 4.b Empty the Receiver Plate and blot with a paper towel.
- 4.c Spin at 300 x g for 3 minutes.

Elute

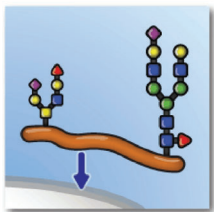
- 5.a Remove the Racks from the Receiver Plates and place on top of a Collection Plate.

NOTE: For a colorimetric measurement (A_{590}/A_{450}) of protein concentration, use the Purification Collection Plate. For direct measurement (A_{280}) of concentration, use a UV-compatible, flat-bottom, half-area plate. If no protein determination will be made, a PCR plate may be used.

- 5.b Pipet 50 μl of Eluent into the Sample Cup of each AssayMAP PA50 Cartridge.
- 5.c Spin at 300 x g for 3 minutes.
- 5.d Remove the Racks and dispose of the Cartridges.

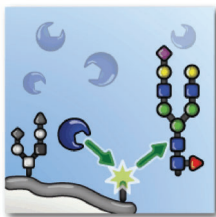
Measure (optional)

6. Measure the absorbance on a plate reader at 280 nm.



Digest

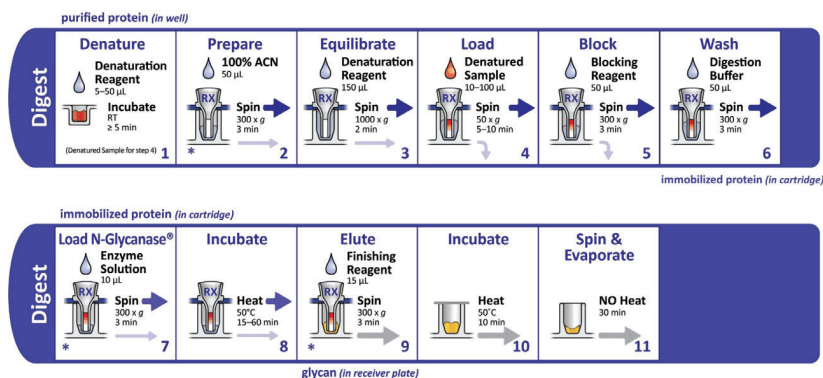
Samples (antibodies or other glycoproteins) are denatured and immobilized.



N-Glycans are released by N-Glycanase and eluted.

Overview

- 1 Denature
- 2 Prepare
- 3 Equilibrate
- 4 Load
- 5 Block
- 6 Wash
- 7 Load N-Glycanase
- 8 Incubate
- 9 Elute (and Finish)
- 10 Incubate
- 11 Spin & Evaporate



Reagents and other Supplies

Glycoprotein Samples

NOTE: The quantity of purified sample loaded to the RX Cartridge may contain NO MORE than 50 μ g total protein, the quantitative binding capacity of the RX Cartridge.

RX Cartridges (supplied with the Kit, 1 per sample)

Prepare two balanced RX Cartridge assemblies (Cartridges on Racks on Receiver Plates with Lids).

Denaturation Reagent (supplied with the Kit)

Acetonitrile (100%, HPLC-grade), 50 μ l/sample

Blocking Reagent (supplied with the Kit)

Digestion Buffer (prepared below)

Enzyme Solution (prepared below)

Finishing Reagent (supplied with the Kit)

Aluminum Sealing Film (supplied with the Kit; film may be cut with scissors if using less than the full Kit)

Preparation of Reagents

Digestion Buffer

NOTE: May be prepared up to one week before use. Store at 2-8°C.

25x Digestion Buffer (supplied with the Kit)

Ultrapure water

Dilute one volume of 25x Digestion Buffer with twenty-four volumes of ultrapure water to obtain Digestion Buffer.

Specifically, add 0.4 ml of 25x Digestion Buffer to 9.6 ml of ultrapure water to make 10 ml of Digestion Buffer.

For fewer samples, prepare 100 μl of Digestion Buffer for each sample to be processed.

Cap tightly and vortex on high for 10 seconds to mix.

Enzyme Solution

NOTE: Should be prepared on the day of use. Store at RT.

N-Glycanase (supplied with the Kit)
Digestion Buffer (prepared above)

Spin the N-Glycanase briefly to collect the contents in the base of the vial, and pipet up and down several times to mix prior to use.

In a separate vial, prepare a mixture of 3 μl of N-Glycanase and 9 μl of Digestion Buffer for each sample to be processed. Pipet up and down several times to mix.

To prepare 24 samples, add 72 μl of N-Glycanase to 216 μl of 1x Digestion Buffer in a pipette basin. Pipet up and down several times to mix.

NOTE: The pipette basin requires a minimum of ~ 100 μl volume, so for fewer than 8 samples, do not use a basin.

Procedure

NOTE: An incubation at elevated temperature is required for full deglycosylation. Before beginning, be sure each Incubation Block has equilibrated to 50°C: a thermometer may be placed in the corner well of Incubation Block to monitor the temperature.

Denature

- 1.a Add Denaturation Reagent to each sample as described in Sample Denaturation (page 7).
- 1.b Pipet up and down to mix.
- 1.c Incubate at room temperature for at least 5 minutes.

NOTE: Proceed through the Prepare, Equilibrate and Load steps without interruption, as evaporation can lead to airlock.

Prepare

- 2.a Pipet 50 μl of 100% acetonitrile into the Sample Cup of each RX Cartridge.
- 2.b Spin at 300 x g for 3 minutes; do not empty the Receiver Plates.

Equilibrate

- 3.a Pipet 150 μl of Denaturation Reagent into the Sample Cup of each RX Cartridge.
- 3.b Spin at 1000 x g for 2 minutes into the same Receiver Plate used for Step 2.b.

NOTE: Do not empty Receiver Plate prior to loading the denatured sample.

Load

4.a Load each Denatured Sample into the Sample Cup of an RX Cartridge (see Sample Loading Technique in the GlykoPrep Guidebook).

4.b Empty the Receiver Plate and blot with a paper towel.

NOTE: Discard waste containing acetonitrile according to waste disposal procedures.

4.c Spin at 50 x g until all Sample Cups are empty. The estimated spin time is 5 minutes for volumes between 10 and 50 μl or 10 minutes for volumes up to 100 μl .

Block

5.a Pipet 50 μl of Blocking Reagent into the Sample Cup of each RX Cartridge.

5.b Empty the Receiver Plate and blot with a paper towel.

5.c Spin at 300 x g for 3 minutes; do not remove the Receiver Plate.

Wash

6.a Pipet 50 μl of Digestion Buffer into the Sample Cup of each RX Cartridge.

6.b Spin at 300 x g for 3 minutes.

Load N-Glycanase

7.a Transfer RX Cartridges to fresh Receiver Plates.

7.b Pipet 10 μl of Enzyme Solution into the Sample Cup of each RX Cartridge.

7.c Spin at 300 x g for 3 minutes;
DO NOT DISCARD FLOW-THROUGH.

Incubate

8. Incubate RX Cartridge assemblies on the equilibrated Incubation Blocks (Heater setting 50°C) for the chosen Incubation Time (not to exceed 60 minutes; see Digestion Module Time, page 8).

Elute (and Finish)

9.a Remove the RX Cartridge assemblies from the Incubation Blocks. Keep the Heater on for the next Incubate step.

NOTE: If condensation is apparent, spin at 300 x g for 3 minutes and tap dish gently on the bench top to release Cartridges that may be stuck to the Lid.

9.b Pipet 15 μl of Finishing Reagent into the Sample Cup of each RX Cartridge; do not remove Rack from Receiver Plate.

9.c Spin at 300 x g for 3 minutes.

9.d Remove Cartridges and Rack from the Receiver Plate. The eluted N-glycans are in the Receiver Plate; DO NOT DISCARD.

Incubate

10. Cover the Receiver Plate with Aluminum Sealing Film and incubate on the equilibrated Incubation Block (with Heater still set at 50°C) for 10 minutes.

NOTE: After removing the Receiver Plate, adjust the temperature setting to 70°C. A thermometer may be placed in the corner thermometer well of the Incubation Block to confirm the temperature.

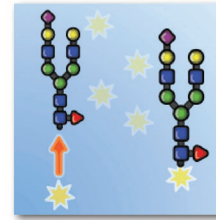
Spin & Evaporate

11. Remove the Aluminum Sealing Film and dry the N-glycans in a centrifugal evaporator (SpeedVac, heat setting turned to the off position) for 30 minutes or until fully dry. This condenses N-glycans into a pellet small enough to be dissolved by 5 μ l of Labeling Reagent in the next step.

NOTE: The Finishing Reagent must be removed promptly by drying. Store N-glycans at -20 or -80°C in the dark if further processing will not take place immediately.

NOTE: Retain the RX Cartridges to recover the deglycosylated protein for further analysis (see Tips & Hints).

NOTE: If N-glycans are to be labeled, now would be a good time to prepare the Labeling Reagent.



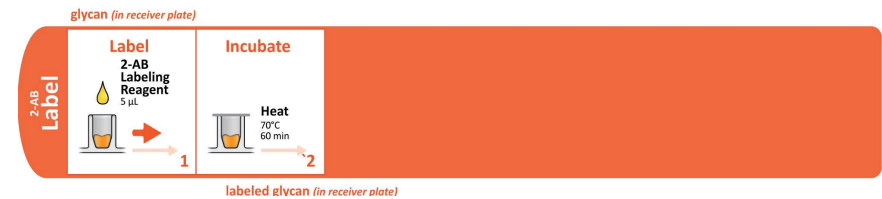
Label

2-AB Labeling Reagent is optimized to label via reductive amination with high labeling efficiency in just one hour.

Overview

Label

Incubate



Reagents and other Supplies

N-Glycan Samples (≤ 30 μ g of dried N-glycans with a free reducing end in a Receiver Plate)

Aluminum Sealing Film (supplied with the Kit; film may be cut with scissors if using less than the full Kit)

2-AB Labeling Reagent (prepared below, no more than one hour before use)

Preparation of Reagents

2-AB Labeling Reagent

NOTE: 2-AB Labeling Reagent should be prepared no more than one hour before use.

NOTE: 2-AB Labeling Reagent components are hazardous. Please refer to the Safety Data Sheets included in the Kit or on our website.

NOTE: Both the 2-AB Solution and Reductant Solution are hygroscopic; minimize exposure to air and protect from exposure to light. The individual reagents may be resealed, repackaged with the desiccant in the resealable bag, and frozen (-20°C) for storage up to 6 months; return to RT before opening for use to minimize condensation.

2-AB Solution (supplied with the Kit)
Reductant Solution (supplied with the Kit)
Aluminum Sealing Film (supplied with the Kit)
Vial, 1-ml, for use with organic solvents

Allow the 2-AB Solution and Reductant Solution vials to come to room temperature in the sealed desiccant bag before removing them. Before opening each vial, flick it or gently tap it on a flat surface to dislodge any liquid adhering to the underside of the cap and ensure that the contents collect at the bottom.

NOTE: Because the viscosity and high surface tension of the 2-AB Labeling Reagent prevent it from spreading out in a pipette basin, do not use a basin for this mixture. If using a multichannel pipette, one column of a clean, empty Receiver Plate or other 96-well plate compatible with organic solvents may be used to hold 2-AB Labeling Reagent, which may then be dispensed into rows of the N-Glycan sample plate.

In a separate vial, prepare a mixture of 3 μ l of 2-AB Solution and 3 μ l of Reductant Solution for each sample to be processed.

Cap tightly and vortex on high for 10 seconds to mix; briefly spin down in a centrifuge.

Tightly cap the 2-AB Solution and Reductant Solution vials, return them to the desiccant-containing bag and store at -20°C.

Procedure

NOTE: Perform this procedure using appropriate personal safety protection, eyeglasses and nitrile gloves.

Label

- 1.a Add 5 μ l of 2-AB Labeling Reagent to the side of each well containing dried N-Glycan Sample. Gently tap the plate on the counter to force the droplet down to the bottom of the well.
- 1.b Seal the plate with Aluminum Sealing Film. Vortex the plate on high for 1 minute to mix and centrifuge (1 minute at 1000 x g) to ensure the samples are collected at the bottom of the wells.

Incubate

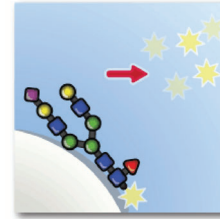
NOTE: Before use, be sure each Incubation Block has equilibrated to 70°C.

- 2.a Incubate sealed plate at 70°C on the equilibrated Incubation Blocks for 1 hour.

NOTE: This is a good time to prepare the 96% Acetonitrile for Cleanup.

- 2.b Remove the Receiver Plate from the Incubation Block and allow it to reach room temperature (5 minutes).
- 2.c In a fume hood, remove the Aluminum Sealing Film. It is normal for condensate to collect on the underside of the film. DO NOT centrifuge the plate to collect the condensate.

Proceed immediately to Cleanup.

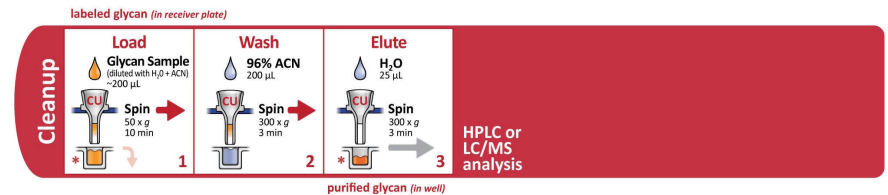


Cleanup

CU Cartridges allow most hydrophobic, non-glycan contaminants to be washed through; N-glycans are then eluted with water.

Overview

- 1 Load
- 2 Wash
- 3 Elute



Reagents and other Supplies

N-Glycan Samples from 2-AB Labeling (≤ 30 µg of N-glycans in 5 µL of 2-AB Labeling Reagent)

CU Cartridges (supplied with the Kit, 1 per sample)
Prepare two balanced CU Cartridge assemblies
(Cartridges on Racks on Waste Plates with Lids)

Acetonitrile (100%, HPLC-grade)

96% Acetonitrile Solution (prepared below)

Ultrapure water

Cleanup Collection Plate (supplied in the AM200 Starter Labware Set, or equivalent)

Preparation of Reagents

96% Acetonitrile Solution

NOTE: May be prepared up to one week before use. Store sealed at room temperature.

Ultrapure water
Acetonitrile (100%, HPLC-grade)

To make 15 ml (enough for a full 24-well Kit) of 96% Acetonitrile Solution (v/v), add 0.6 ml of ultrapure water to a glass, graduated cylinder. Bring the volume up to 15 ml with 100% acetonitrile. Transfer to a glass storage vessel, cap tightly and swirl gently to mix.

Scale as necessary (prepare 250 μ l per sample).

Procedure

NOTE: DO NOT use Receiver Plates in this procedure; build the stack with Waste Plates (~450- μ l well volume) instead. This entire section is performed with the CU Cartridges “tips free.”

Load

- 1.a Add 20 μ l of ultrapure water to each N-Glycan Sample.
- 1.b Add 180 μ l of 100% Acetonitrile to each N-Glycan Sample. Pipet up and down to mix.
- 1.c Transfer each N-Glycan Sample into the Sample Cup of a CU Cartridge. This must be done quickly because Acetonitrile has very low viscosity and may drip from the pipette tip; each sample may be pipetted in multiple rounds in order to achieve a quantitative transfer.

NOTE: Air bubbles are not a concern with this concentration of Acetonitrile.

1.d Spin at 50 x g for 10 minutes or until the CU Cartridges are empty.

1.e Empty the Waste Plate.

NOTE: Discard waste containing acetonitrile and 2-AB Labeling Reagent according to waste disposal procedures.

Wash

2.a Pipet 200 μ l of 96% Acetonitrile Solution into the Sample Cup of each CU Cartridge.

2.b Spin at 300 x g for 3 minutes.

2.c Empty the Waste Plate.

Elute

3.a Place each racked set of CU Cartridges over a clean Cleanup Collection Plate.

NOTE: Because the eluate contains traces of organic solvent, polystyrene plates should NOT be used. Any polypropylene ANSI/SBS 96-well microplate may be used as a collection plate. To facilitate complete product recovery, we recommend plates with conical bottoms, such as PCR plates or the Cleanup Collection Plates provided in the AM200 Starter Labware Set.

3.b Pipet 25 μ l of ultrapure water into the Sample Cup of each CU Cartridge.

NOTE: Up to 200 μ l of water may be used if more dilute N-glycans are desired.

NOTE: To protect N-glycans for long-term storage, an aqueous buffer compatible with the intended analysis method may be used instead of water.

3.c Spin at 300 x g for 3 minutes.

The Cleanup Collection Plate now contains the purified N-glycans; DO NOT DISCARD.

Mix the eluate by pipette action or vortexing prior to analysis to ensure homogeneity. N-Glycan Samples are now ready to be analyzed. If not analyzed immediately, store sealed at -20°C in the dark.

ANALYSIS OF LABELED N-GLYCANS

Use standard techniques, such as Liquid Chromatography (LC), Mass Spectrometry (MS), or a combination of the two, to analyze the aqueous eluate containing eluted, labeled N-glycans (see Tips and Hints, below).

TIPS & HINTS

Optimizing Excitation/Emission Wavelengths

Optimal excitation/emission wavelengths for 2-AB Dye conjugated to an N-glycan may vary depending upon the optical configuration of the instrument used.

Excitation/emission pairs that have been used together include 250/428 nm (Melmer et al., 2010), 330/420 nm (Bigge et al., 1995) and 360/428 nm (used by ProZyme with the Waters Acquity UPLC®; Haxo et al., 2012).

Recovery of the Deglycosylated Protein from the Digestion (RX) Cartridge

Often, the deglycosylated protein is analyzed to evaluate the completeness of deglycosylation using such electrophoretic methods as SDS-PAGE or microfluidic lab-on-a-chip technology. Please contact us for guidelines for eluting your glycoprotein from the RX Cartridge.

Calculating the Mass of Glycans Labeled with 2-AB

The reductive amination reaction results in the loss of an oxygen atom. The mass of the 2-AB-labeled N-glycan is obtained using the following formula:

$$\text{Mass}_{\text{Glycan}} + \text{Mass}_{2\text{-AB}} = \text{Mass}_{2\text{-AB-Labeled Glycan}} - \text{Mass}_{\text{Oxygen}}$$

Mass Added to Glycan

Monoisotopic 120.06875

Average 120.2

Direct LC Analysis of N-Glycans After Elution

If N-Glycan Samples will be analyzed by LC directly following elution from the CU Cartridges, the 96-well, polypropylene plate with a pierceable lid available from MicroLiter Analytical Supplies (cat# 07-1211N) may serve as a Collection Plate.

Alternatively, Cleanup Collection Plates may be heat sealed with pierceable foil (e.g., Thermo Easy Pierce 20 μm Foil, #AB-1720) using a microplate heat sealer (e.g., Thermo ALPS 50 V Semi automated Microplate Heat Sealer, #AB-1443).

REFERENCES

- Bigge, J. C., T. P. Patel, J. A. Bruce, P. N. Goulding, S. M. Charles, R. B. Parekh. Nonselective and Efficient Fluorescent Labeling of Glycans Using 2-Amino Benzamide and Anthranilic Acid. *Anal Biochem* 230: 229–38 (1995).
- Melmer, M., T. Stangler, M. Schiefermeier, W. Brunner, H. Toll, A. Ruppachter, W. Lindner and A. Premstaller. HILIC Analysis of Fluorescence-Labeled N-Glycans from Recombinant Biopharmaceuticals. *Anal Bioanal Chem* 398: 905–914 (2010).
- Haxo, T., J. Hyche, M. Kimzey, S. Lockhart, C. Nishida, S. Pourkaveh, J. Wegstein, Y. Q. Yu and D. J. Phillips. An Integrated Strategy for N-Glycan Sample Preparation and Analysis Suitable for All Stages of Therapeutic Protein Discovery, Characterization, Manufacture and Quality Release. Poster session presented at: Well Characterized Biotechnology Pharmaceuticals 16th Symposium on the Interface of Regulatory and Analytical Sciences for Biotechnology Health Products 2012 Jan 22–25 San Francisco, CA, USA.

Visit ProZyme's website for additional information, downloadable posters and instructional videos:

<http://www.prozyme.com/glykoprep>

TechNote TNGP100 GlykoPrep Guidebook - General tips, tricks and troubleshooting suggestions when using Kits or modules:

<http://www.prozyme.com/documents/TNGP100.pdf>

TECHNICAL ASSISTANCE

ProZyme is committed to developing rapid, automatable methods for glycan analysis. Call us to discuss products in development.

If you have any questions or experience difficulties regarding any aspect of our products, please contact us:

TOLL FREE **(800) 457-9444** (US & CANADA)
PHONE **(510) 638-6900**
FAX **(510) 638-6919**
E-MAIL **info@prozyme.com**
WEB **www.prozyme.com**

ProZyme values customer opinions and considers customers an important source for information regarding advanced or specialized uses of our products. We encourage you to contact us. We welcome your suggestions about product performance or new applications and techniques.

LICENSE TO USE

Purchase of this ProZyme product (the "Product") from ProZyme or one of its authorized distributors comes with a limited, non-transferable Use License to the purchaser ("Purchaser") under ProZyme patents (the "ProZyme Patents"), which Use License is subject to the following terms and conditions.

1. The Use License is only for Purchaser's in-house, in-vitro use of the purchased Product. Purchaser is not licensed to use, and shall not use the Product in any in vivo application and, for avoidance of doubt, is specifically prohibited from any use of the Product in humans.
2. The Product shall not be made available by Purchaser to any third party in any form, separately or in combinations, whether for any monetary or other consideration or at no charge, with the following exception: the Product may be made available to a third party who (a) agrees to be bound by all the terms and restrictions of this Agreement and (b) agrees to use it only for purposes of evaluation.
3. This Use License does not include the right to use the Product to perform for, or offer commercial services of any kind to, a third party, such as, without limitation, reporting the results of analyses or other activities to the third party for a fee or other commercial consideration. Persons wishing to perform such services must obtain a separate Commercial Services License from ProZyme, Inc.
4. Purchaser is responsible for qualification of the Product for the Purchaser's specific use.

Certain ProZyme products include components or reagents obtained under license from other companies. Purchase of these ProZyme products from ProZyme or one of its authorized distributors includes a license from the originating company to use the component(s) or reagent(s), but only (1) with the other materials included in the ProZyme product, and (2) for the purposes specified on ProZyme's product, on its packaging, or in the accompanying literature.

For a complete list of ProZyme Patents, please visit our website at

<http://www.prozyme.com/patent.html>

The AssayMAP technology is licensed from Agilent Technologies.

TRADEMARKS AND TRADENAMES

ProZyme, GlykoPrep, Rapid-Reductive-Amination and N-Glycanase are trademarks or registered trademarks of ProZyme, Inc. in the United States and other countries.

Diamond is a registered trademark of Gilson Inc., Middleton, WI, USA.

SpeedVac is a registered trademark of Thermo Fisher Scientific Inc.

Milli-Q is a registered trademark of Millipore Corporation in the United States and/or other countries.

Waters, Acquity and UPLC are registered trademarks of Waters Corporation, Milford, MA, USA.

PRODUCT USE AND WARRANTY

Terms and conditions of sale may be found at:

<http://www.prozyme.com/terms.html>

OTHER PROZYME PRODUCTS & KITS

A wide variety of glycoanalysis products are available from ProZyme. A complete listing is accessible on our website by clicking on GlykoPrep® Rapid Sample Preparation Platform:

<http://www.prozyme.com>

ORDERING INFORMATION

For North American destinations: telephone orders may be placed between 8:00 am and 5:00 pm Pacific Time. Telefax or e-mail orders may be sent or messages recorded anytime.

TOLL FREE (800) 457-9444 (US & CANADA)

PHONE (510) 638-6900

FAX (510) 638-6919

E-MAIL info@prozyme.com

WEB www.prozyme.com

Outside North America:

A list of ProZyme's distributors, with contact information, may be found at:

<http://www.prozyme.com/distributors.html>

If there is no distributor in your area, instructions for placing an international order may be found at:

<http://www.prozyme.com/ordering.html>



3832 Bay Center Place
Hayward, CA 94545-3619
USA

TOLL FREE (800) 457-9444

PHONE (510) 638-6900

FAX (510) 638-6919

E-MAIL info@prozyme.com

WEB www.prozyme.com

Rev. AD