

Agilent Buffer Advisor

User's Guide



Agilent Technologies

Notices

© Agilent Technologies, Inc. 2012

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.

Manual Part Number

G5617-90000

Edition

5/12

Printed in Germany

Agilent Technologies
Hewlett-Packard-Strasse 8
76337 Waldbronn

This product may be used as a component of an in vitro diagnostic system if the system is registered with the appropriate authorities and complies with the relevant regulations. Otherwise, it is intended only for general laboratory use.

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

If software is for use in the performance of a U.S. Government prime contract or subcontract, Software is delivered and licensed as “Commercial computer software” as defined in DFAR 252.227-7014 (June 1995), or as a “commercial item” as defined in FAR 2.101(a) or as “Restricted computer software” as defined in FAR 52.227-19 (June 1987) or any equivalent agency regulation or contract clause. Use, duplication or disclosure of Software is subject to Agilent Technologies’ standard commercial license terms, and non-DOD Departments and Agencies of the U.S. Government will

receive no greater than Restricted Rights as defined in FAR 52.227-19(c)(1-2) (June 1987). U.S. Government users will receive no greater than Limited Rights as defined in FAR 52.227-14 (June 1987) or DFAR 252.227-7015 (b)(2) (November 1995), as applicable in any technical data.

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

1	Introduction to Buffer Advisor	5
	What is Buffer Advisor?	6
2	A Short Excursion into Ion Exchange Chromatography	9
	Properties of Proteins	10
	Ion Exchange: Salt Gradient	12
	Ion Exchange: pH Gradient (Chromatofocusing)	14
3	Columns and Buffers	17
	Column Selection	18
	Buffer Selection	20
4	Starting with Buffer Advisor	23
	Installation	24
	User Interface Overview	25
5	Ion Exchange Experiments	31
	Salt Gradient with Single Buffer System	32
	Short pH Gradient with Single Buffer System	34
	pH Gradient with Multi-Buffer System	36
	pH Scouting with Single Buffer System	38
	Generate and export a gradient time table	40
	Import a gradient timetable into the G5611A pump driver	41
6	Troubleshooting and Tips	43
	Messages shown by buffer advisor	44
	Essential measurement practices	46
7	Validation Tests and Specifications	49

Contents



1

Introduction to Buffer Advisor

What is Buffer Advisor? 6

This chapter provides an overview of the Agilent Buffer Advisor software.



What is Buffer Advisor?

The Agilent Buffer Advisor Software is a calculator that helps you to set up a salt gradient or pH gradient for ion exchange chromatography. In general, any aqueous buffer can be generated from up to 4 stock solutions. This also includes optimization of ionic strength, or pH scouting for SEC, HIC and Affinity Chromatography

The use of Buffer Advisor typically involves the following steps:

- Definition of buffer concentration, pH and maximum salt concentration for an ion exchange separation, and selection of the most suitable concentrations of stock solutions.
- Verification that the selected buffer system is suitable for the separation task.
- Automatic generation of a corrected pump timetable for a salt gradient or pH gradient.

The mixing principle of Agilent Buffer Advisor is depicted in [Figure 1](#) on page 6 for salt gradients, and [Figure 2](#) on page 7 for pH gradients.

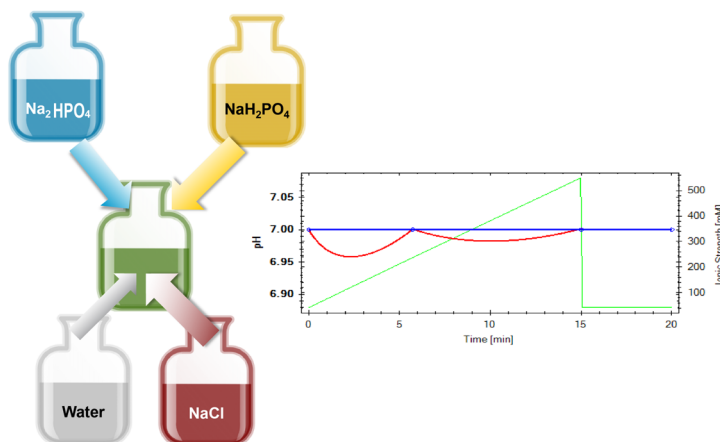


Figure 1 Quaternary mixing of four stock solutions (water, salt solution, acid and base) with the Agilent Bio-inert LC and Buffer Advisor generates a linear salt gradient at constant pH.

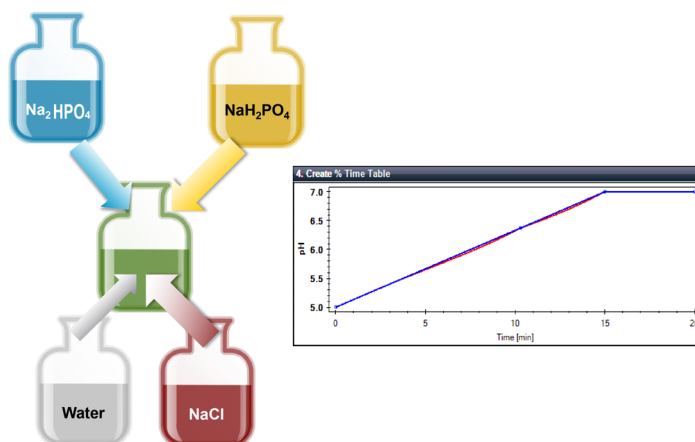


Figure 2 Proportioning of three solutions (water, acid, base) with Agilent Bio-inert quaternary LC and Buffer Advisor results in formation of a linear pH gradient.

Quaternary solvent mixing using the Agilent Buffer Advisor provides the capability of automatically forming several mobile phases of different pH values for different methods from four stock solvents. This enables automated method development (pH scouting) by screening mobile phases for an analyte separation at different pH values without the necessity to prepare the different buffers manually. In addition, the Agilent Buffer Advisor software corrects a gradient timetable with compensation points (added timetable steps), such that the pH is kept constant while the ionic strength is changing.

The corrected timetable can be imported into the Agilent Bio-inert pump (G5611A) driver and is displayed as a pump timetable in the user interface of the respective ChemStation or EZChrom method.

1 Introduction to Buffer Advisor

What is Buffer Advisor?



2

A Short Excursion into Ion Exchange Chromatography

Properties of Proteins 10

Ion Exchange: Salt Gradient 12

Ion Exchange: pH Gradient (Chromatofocusing) 14

This chapter provides a short introduction to the challenges posed by ion exchange chromatography and how you can use the Agilent Buffer Advisor software to solve them.



Properties of Proteins

Proteins consist of amino acids with both acidic and basic functional groups, leading to a net positive, neutral or negative surface charge. Because of differences in pK_a of the basic and acidic amino acids, the net charge of the molecule is dependent on the pH of the surrounding medium. At high pH, the net charge of the protein changes towards negative, whereas at low pH, it becomes positive. The pH where the net charge of the protein is zero is called the *isoelectric point*, pI .

In *ion exchange chromatography*, two approaches can be used to exploit the net charge differences of proteins for separation:

- Most commonly used are *salt gradients*, where increasing ionic strength leads to the separation of proteins.
- Alternatively, *pH gradients* can be used to avoid the necessity of applying highly concentrated salt solutions.

In ion exchange chromatography, knowledge of a protein's pI is used to select the starting conditions to ensure that the protein interacts with the stationary phase. A mobile phase pH below the pI of the protein of interest allows the use of a Cation Exchanger column, whereas a mobile phase pH above the pI enables protein separation on an Anion Exchanger column (Figure 3 on page 11). In both cases, the protein binds to the column according to its current charge. A negatively charged column is used in Cation Exchange, whereas a positively charged column is used in Anion Exchange.

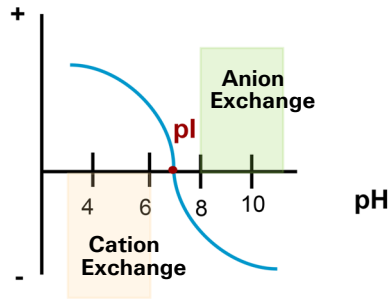


Figure 3 Isoelectric point (pI) of a theoretical protein of 7. If the protein is stable above its pI, Anion Exchange is recommended, if it is stable below its pI, Cation Exchange is recommended.

When a *salt gradient* is used, the ionic strength of the eluent is increased to elute the proteins in the order of increasing binding strength of the column.

If a *pH gradient* is used, the proteins elute at the pH where the net charge is zero, at their isoelectric points.

[Table 1](#) on page 11 shows a list of proteins and their pIs. A buffer system and an appropriate column are selected for the analytical task depending on the pI of the protein. As a rule of thumb, proteins with high pI are separated on cation-exchange columns, proteins with low pI on anion exchanger columns.

Table 1 List of example proteins and their pIs

Protein	pI	Source
Amyloglucosidase	3.6	Aspergillus niger
Ovalbumin	4.7	Chicken, egg white
BSA	4.9	
Human insulin	5.4	
Bovine carbonic anhydrase	6.5, 5	Cow, erythrocyte
Myoglobin	7.4, 6.9	Horse, muscle
Ribonuclease A	8.88	Cow, pancreas
Cytochrome C	10.7, 9	Horse, heart
Lysozyme	11.3	Chicken, egg white

Ion Exchange: Salt Gradient

A salt gradient triggers a *competing* process. In cation exchange chromatography, positively charged proteins bind to a negatively charged solid support. The mobile phase buffer is kept at a constant pH below the protein's pI to keep the protein's net charge positive. Salt ions (for example, NaCl) are used as competing agents so that at a low concentration of salt (low ionic strength) the protein binds to the column, whereas with increasing salt concentration the positively charged sodium ions displace the positively charged proteins, leading to elution. The higher the net charge of the protein, the later in the gradient it will elute.

Binding

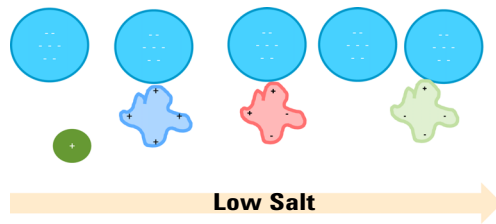


Figure 4 In a first step, positively charged proteins are bound to the column.

Elution

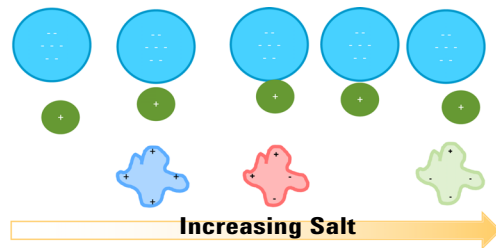


Figure 5 The salt concentration is increased and proteins elute.

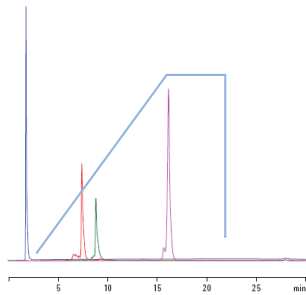


Figure 6 Example chromatogram of Cation Exchange chromatography employing a salt gradient. Ovalbumin, Ribonuclease A, Cytochrome c, Lysozyme were separated in pH 6.5 phosphate buffer. Protein separation with a salt gradient from 0–800 mM NaCl, 20 mM phosphate buffer, pH 6.5, automatically generated by the Agilent Buffer Advisor software.

Table 2 Conditions

Columns	Agilent Bio WCX NP10, 4.6 × 250 mm SS Agilent Bio WCX NP5, 4.6 × 250 mm SS
Mobile phase	A: Water B: 1.6 M NaCl C: 40.0 mM NaH ₂ PO ₄ D: 40.0 mM Na ₂ HPO ₄ By combining predetermined proportions of C and D, 20 mM buffer solutions at the desired pH range were produced.
Gradient	0 – 50 % B, 0 – 20 min 50 % B, 20 – 25 min 0 % B, 25 – 35 min
Flow rate	1.0 mL/min
Temperature	Ambient
Sample	Ovalbumin, Ribonuclease A, Cytochrome c, Lysozyme
Detection	UV, 220 nm
Instrument	Agilent 1260 Infinity Bio-inert HPLC system

Ion Exchange: pH Gradient (Chromatofocusing)

A pH gradient consists of a starting buffer and an elution buffer. In the starting buffer, the net surface charge of the protein should be such that it can interact with the functional group of the ion exchanger resin. For example, in cation exchange, the starting pH of the buffer system is low. The pH is then increased linearly, resulting in de-protonation of the protein's amino acids, and subsequently to elution once the pI of the protein and the pH of the mobile phase match. At the same time, the buffer forms a pH gradient across the column. Proteins with less positive charges elute earlier than those with more positive net charges (see [Figure 7](#) on page 14).

In contrast, in anion exchange chromatography, the starting conditions are at high pH, and the pH decreases towards acidic during the chromatographic run.

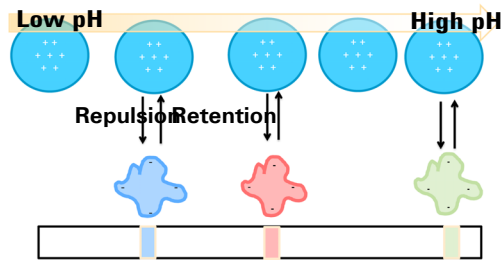


Figure 7 Example of a pH gradient: Proteins with same pI are focused in bands on the column.

The advantages of IEX with a pH gradient are threefold:

- improved resolution
- the pI of the protein can be correlated to the pH of the buffer system
- salt solutions can be avoided (reducing ionic interactions and preventing corrosion)

A pH gradient is preferable, but it is difficult to form a linear pH gradient over a wide pH range because a single buffer system can provide sufficient buffer capacity over only a narrow pH range.

The Buffer Advisor software is an ideal tool that helps to select a suitable buffer system for a specific separation task.

2 A Short Excursion into Ion Exchange Chromatography

Ion Exchange: pH Gradient (Chromatofocusing)



3

Columns and Buffers

Column Selection 18

Buffer Selection 20

This chapter gives some background information on the selection of columns and buffers for ion exchange chromatography.



Column Selection

The column and separation buffer are selected depending on the charge of the protein. In ion exchange chromatography, the pH of the mobile phase buffer must be between the pI or pKa of the charged molecule and the pKa of the charged groups on the stationary phase.

In *anion exchange chromatography*, a strong or a weak anion exchange column containing a quaternary ammonium ion, or a weak anion exchanger with either a tertiary or secondary amine functional group such as DEAE (diethylaminoethyl) is used.

cation exchange chromatography is conducted with either a strong or a weak cation exchange column containing a sulfonium ion, or with a weak cation exchanger usually having a carboxymethyl (CM) functional group. A counter-ion, often Na⁺ maintains electroneutrality.

Agilent offers a wide range of columns for ion exchange:

Table 3 Selection of high resolution ion-exchange columns for charged-based protein separation

Column name	Dimensions	Particle size (µm)	Product number
Agilent Bio MAb	4.6 × 50 mm	3	5190-2403
Agilent Bio MAb	4.6 × 250 mm	5	5190-2407
Agilent Bio MAb	4.6 × 50 mm	5	5190-2408
Agilent Bio MAb	2.1 × 250 mm	5	5190-2411
Agilent Bio MAb	2.1 × 50 mm	5	5190-2412
Agilent Bio SCX	4.6 × 50 mm	3	5190-2423
Agilent Bio SCX	4.6 × 250 mm	5	5190-2427
Agilent Bio SCX	4.6 × 50 mm	5	5190-2428
Agilent Bio WCX	4.6 × 50 mm	3	5190-2443
Agilent Bio WCX	4.6 × 250 mm	5	5190-2447
Agilent Bio WCX	4.6 × 50 mm	5	5190-2448

Table 3 Selection of high resolution ion-exchange columns for charged-based protein separation

Column name	Dimensions	Particle size (μm)	Product number
Agilent Bio SAX	4.6 \times 50 mm	3	5190-2463
Agilent Bio SAX	4.6 \times 250 mm	5	5190-2467
Agilent Bio SAX	4.6 \times 50 mm	5	5190-2468
Agilent Bio WAX	4.6 \times 50 mm	3	5190-2483
Agilent Bio WAX	4.6 \times 250 mm	5	5190-2487
Agilent Bio WAX	4.6 \times 50 mm	5	5190-2488

More information can be found at
<http://www.chem.agilent.com/en-US/Products/columns-supplies/lc-lc-mscolumns>.

Buffer Selection

Typically, low ionic strength buffers are used in ion exchange chromatography so that the charged proteins are retained by the stationary phase that has the opposite charge. Proteins with the same charge as the stationary phase will simply flow through without binding. The ion exchange matrix is washed with additional low ionic strength buffer to completely wash out any remaining unbound species, and the bound species are differentially eluted by buffers containing increasing concentration of salt. As the ionic strength of the mobile phase increases, salt ions compete for binding to the charges on the ion exchange matrix, displacing the bound macromolecule and allowing them to elute from the matrix.

Therefore anionic (negatively-charged) buffers are used for *cation exchange* and cationic (positively-charged) buffers are used for *anion exchange*.

A buffer is characterized by its pKa, the negative logarithm of the acid dissociation constant (Ka). Buffer capacity is highest when the pH equals the pKa of the buffer.

- If the pH of the buffer stays below the pI then the net protein charge is positive.
- If the pH of the buffer stays above the pI then the net protein charge is negative.
- With salt gradient elution, start at least one pH unit away from the protein's pI.
- Use a buffer concentration that is sufficient to maintain buffer capacity to keep pH constant, usually between 20 and 50 mM.

Table 4 Buffers for Cation Exchange Chromatography

Name	pKa (25 °C)
Citric Acid	3.13
Lactic Acid	3.86
Succinic Acid	4.21
Acetic Acid	4.75

Table 4 Buffers for Cation Exchange Chromatography

Name	pKa (25 °C)
Methyl Malonic Acid	5.76
MES	6.27
Phosphate	7.20
HEPES	7.56
BICINE	8.33

Table 5 Buffers for Anion Exchange Chromatography

Name	pKa (25 °C)
N-Methyl-Piperazine	4.75
Piperazine	5.33
Bis-Tris	6.48
Bis-Trispropane	6.65 9.1
Triethanolamine	7.76
TRIS	8.07
N-Methyldiethanolamine	8.53
Propane 1,3 diamino	8.88
Ethanolamine	9.50
Piperazine	9.73
Propane 1,3 diamino	10.55
Piperidine	11.12

3 Columns and Buffers

Buffer Selection



4

Starting with Buffer Advisor

Installation	24
User Interface Overview	25
Buffer Selection Area	26
List of Available Buffers	27
Recipes	28

This chapter tells you how to install the Agilent Buffer Advisor, and gives you an introduction to the different parts of the graphical user interface.



Installation

- 1 Place the CD-ROM in your CD/DVD drive.
- 2 Double-click the .exe file.

Buffer Advisor will be automatically installed in the Windows default program folder.

To verify that the installation was successful, go to the respective folder, right-click on the file, select **Properties** and verify that the version A.01.01 is correct.

The software starts in demo mode. You can use all capabilities of the Buffer Advisor software but you cannot create a gradient result timetable and save it. To use all the features of Buffer Advisor, enter the license key number under **Help**.

User Interface Overview

The screenshot displays the Agilent Lab Advisor software interface, divided into four main sections:

- 1. Select Buffer & Gradient Mode:** This section allows users to choose between 'Single Buffer (pH / Salt Gradient)' and 'Composite Buffer (Wide Range pH Gradient)'. It also includes options for 'Cation Exchange' and 'Anion Exchange', and a dropdown menu for 'Sodium Phosphate (NaH2PO4+Na2HPO4) - 7.2'.
- 2. Define Gradient Table:** A table with columns for Time, Salt, pH, and Buffer. The data shown is:

Time	Salt	pH	Buffer
0	0	7	20
15	500	7	20
15.01	0	20	20
20	0	20	20
- 3. Compose Stock Solutions:** This section lists stock solutions A, B, C, and D with their respective concentrations and recommended values.

Solution	Concentration	Recommended
A: Water	1500.0 mM	1500.0 mM
B: NaCl	25.0 mM	25.0 mM
C: NaH2PO4	41.0 mM	25.0 mM
D: Na2HPO4	41.0 mM	41.0 mM
- 4. Create % Time Table:** A graph showing pH and Ionic Strength over time. The x-axis is Time [min] (0 to 20), and the y-axis is pH (6.90 to 7.05) and Ionic Strength [mM] (0 to 500). The graph shows a step change in pH and a linear increase in Ionic Strength. A table below the graph shows the resulting gradient data.

Result Pump Gradient Time Table

Time	% A	% B	% C	% D	Init. pH	Calc. pH	IS	BC	Cond.	Status
0	36.6	0	37.5	25.9	7	7	41.2	12	0.252	Correct
5.73	28.6	12.7	25.5	33.2	7	7	238	10.1	2.15	Correct
15	10.8	33.3	18.2	37.7	7	7	550	8.15	4.95	Correct
15.01	36.6	0	37.5	25.9	7	7	41.2	12	0.252	Correct
20	36.6	0	37.5	25.9	7	7	41.2	12	0.252	Correct

Messages

Minimum Buffering Capacity (BC) = 8.15 mM

Area 1: Select Buffer and Gradient Solution

Choose between *single buffer mode*, usually applied for salt gradients, and *composite buffer mode* for pH gradients, and whether you want to perform Anion- or Cation-exchange chromatography.

Area 2: Enter Gradient Timetable

Define the gradient parameters such as time, maximum salt concentration (for salt gradients), pH and buffer concentration.

Area 3: Review Composition of Stock Solutions and Adjust

Concentrations for stock solutions are shown, and the recipe button shows amounts of chemical compound to be added to water to prepare the respective stock solutions. You can also enter your own concentrations of stock solutions and the resulting gradient timetable is calculated accordingly.

Area 4: Gradient Display Section: Process Gradient and Review Optimized Gradient with Compensation Steps

When data is entered, the software calculates the percentages of stock solutions for each channel of the quaternary pump. Furthermore, it calculates whether the pH, salt concentration and buffer concentration entered is suitable for the buffer system that was selected. The timetable shows also additional data, such as buffering capacity or ionic strength. A message area gives hints on optimizing the gradient entered in Area 2.

Buffer Selection Area

1. Select Buffer & Gradient Mode

New... Open... Save...

Single Buffer (pH / Salt Gradient) Composite Buffer (Wide Range pH Gradient)

Cation Exchange Anion Exchange

Sodium Phosphate (NaH₂PO₄+Na₂HPO₄)

In the buffer selection area you can select the type of analysis:

- **Single buffer** is a buffer consisting of an acidic and basic component. For salt gradients, NaCl can be added in increasing concentrations.
- **Composite buffer** is a buffer consisting of multiple buffering components. The composite buffer can be used over a wide pH range without compromising the buffer capacity.

Selection of **Cation/Anion exchange** leads to a dropdown list of buffers for this application. For all buffers, the pKa is displayed, facilitating selection of a buffer related to the analyte's pI, usually near ± 0.5 of a buffer's pKa.

Clicking **New** shows all available buffers in the Agilent Buffer Advisor software, sorted by the above criteria and pKa.

List of Available Buffers

Single buffers

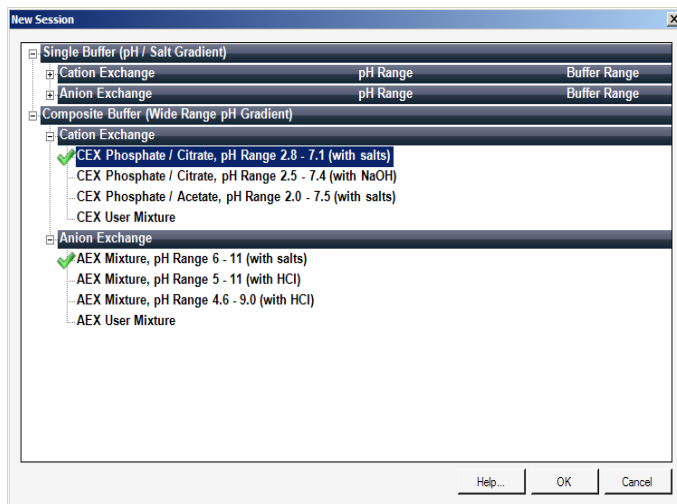
Buffer Name	pH Range	Buffer Range
<input checked="" type="checkbox"/> Sodium Phosphate (NaH ₂ PO ₄ +Na ₂ HPO ₄)	6.1-7.2	7.5-125
<input type="checkbox"/> Sodium Phosphate (H ₃ PO ₄ +NaOH)	2.6-3.6, 5.9-7.4, 10.6-11.1	10-15
<input type="checkbox"/> Sodium Phosphate (H ₃ PO ₄ +Na ₂ HPO ₄)	2.5-3.4, 5.9-7.1	7.5-125
<input type="checkbox"/> Sodium Phosphate (H ₃ PO ₄ +Na ₃ PO ₄)	2.6-3.6, 6.0-7.4, 10.5-11.0	7.5-125
<input checked="" type="checkbox"/> Sodium Citrate (Citric + Tri-Sodium Citrate)	3.0-5.5, 5.5-6.0	7.5-125
<input type="checkbox"/> Sodium Citrate (Citric + NaOH)	2.9-3.7, 3.7-6.2	7.5-15
<input checked="" type="checkbox"/> Formic/Na (acid + Na salt)	3.2-4.4	7.5-125
<input type="checkbox"/> Formic/Na (acid + NaOH)	3.3-4.6	10-50
<input checked="" type="checkbox"/> Lactic/Na (acid + Na salt)	3.2-4.5	7.5-125
<input type="checkbox"/> Lactic/Na (acid + NaOH)	3.4-4.7	7.5-50
<input checked="" type="checkbox"/> Acetic/Na (Acetic+Acetate/Na)	3.9-5.4	7.5-125
<input type="checkbox"/> Acetic/Na (Acetic+NaOH)	4.1-5.6	7.5-50
<input checked="" type="checkbox"/> Succinic/Na (acid + Na salt)	3.6-5.6	7.5-125
<input type="checkbox"/> Succinic/Na (acid + NaOH)	3.9-6.3	10-20
<input checked="" type="checkbox"/> Malonic/Na (acid + Na salt)	2.8-5.5	7.5-125
<input type="checkbox"/> Malonic/Na (acid + NaOH)	2.9-5.5	7.5-25
<input checked="" type="checkbox"/> MES/Na (MES+MES/Na)	5.2-7.1	7.5-125
<input type="checkbox"/> MES/Na (MES+NaOH)	5.5-7.3	7.5-40

Composite Buffers

Six composite buffer mixtures have been predefined:

4 Starting with Buffer Advisor

User Interface Overview



Buffers marked with a green check mark are validated, see “[Validation Tests and Specifications](#)” on page 49.

Depending on the buffer, you select the stock solutions that are displayed on the right-hand side (area 3)

Recipes

Recipe provides the amount of chemicals needed to be weighed in to prepare the stock solutions:

2. Define Gradient Table

Time	Salt	pH	Buffer
0	0	7	20
15	500	7	20
15.01	0	7	20
▶ 20	0	7	20
*			

3. Compose Stock Solutions

A: Water		Recommended
B: NaCl	<input type="text" value="1700"/>	1700 mM
C: NaH2PO4	<input type="text" value="23.5"/>	23.5 mM
D: Na2HPO4	<input type="text" value="38.5"/>	38.5 mM
	<input type="button" value="Recipe..."/>	<input type="button" value="Set"/>

Stock Solution Recipes

Bottle B	<input type="text" value="NaCl: Sodium chloride"/>	Weight 99.348 g and fill up to 1 L.
Bottle C	<input type="text" value="NaH2PO4: Monosodium phosphate"/>	Weight 2.8195 g and fill up to 1 L.
Bottle D	<input type="text" value="Na2HPO4: Sodium phosphate dibasic"/>	Weight 5.4655 g and fill up to 1 L.

4 Starting with Buffer Advisor

User Interface Overview



5 Ion Exchange Experiments

Salt Gradient with Single Buffer System	32
Short pH Gradient with Single Buffer System	34
pH Gradient with Multi-Buffer System	36
pH Scouting with Single Buffer System	38
Generate and export a gradient time table	40
Import a gradient timetable into the G5611A pump driver	41

This chapter describes detailed workflows for four typical ion exchange experiments.



Salt Gradient with Single Buffer System

A salt gradient experiment is usually designed with a buffer around the pKa of the protein of interest. In this example, phosphate buffer (pKa 7.2) is used with a NaCl salt gradient (final concentration 500 mM).

- 1 In the **Select Buffer and Gradient Mode** panel, select the **Single Buffer (pH/Salt Gradient)** option.
- 2 Select the **Cation Exchange** option.
- 3 Open the dropdown menu and select **Sodium Phosphate (NaH₂PO₄+Na₂HPO₄)** from the list of available buffers.
- 4 In the **Define Gradient Timetable** panel, enter the following gradient timetable parameters:

Time (min)	Salt	pH	Buffer
0	0	7	20
15	500	7	20
15.01	0	7	20
20	0	7	20

This timetable keeps the pH close to the pKa, and creates a salt gradient while keeping the pH constant.

- 5 To find the best concentration of stock solutions for this experiment, click **Set** under the heading **Recommended** at the right of the **Compose Stock Solutions** panel.

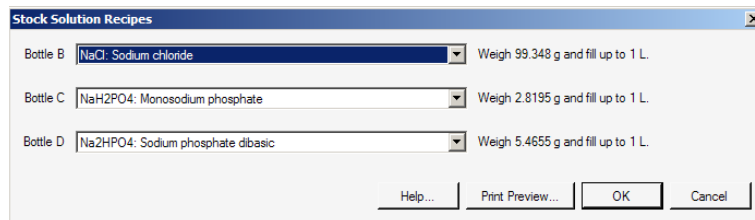
The recommended values above are entered into the fields at the left of the panel.

The screenshot shows three panels from a software interface:

- 1. Select Buffer & Gradient Mode:** Includes buttons for 'New...', 'Open...', and 'Save...'. It has radio buttons for 'Single Buffer (pH / Salt Gradient)' (selected) and 'Composite Buffer (Wide Range pH Gradient)'. Below are radio buttons for 'Cation Exchange' (selected) and 'Anion Exchange'. A dropdown menu shows 'Sodium Phosphate (NaH₂PO₄+Na₂HPO₄)' selected.
- 2. Define Gradient Table:** A table with columns 'Time', 'Salt', 'pH', and 'Buffer'. It contains the same data as the table above.
- 3. Compose Stock Solutions:** Lists components A: Water, B: NaCl, C: NaH₂PO₄, and D: Na₂HPO₄. To the right, under 'Recommended', are input fields with values 1700, 23.5, and 38.5, and a 'Recipe...' button. A 'Set' button is at the bottom right.

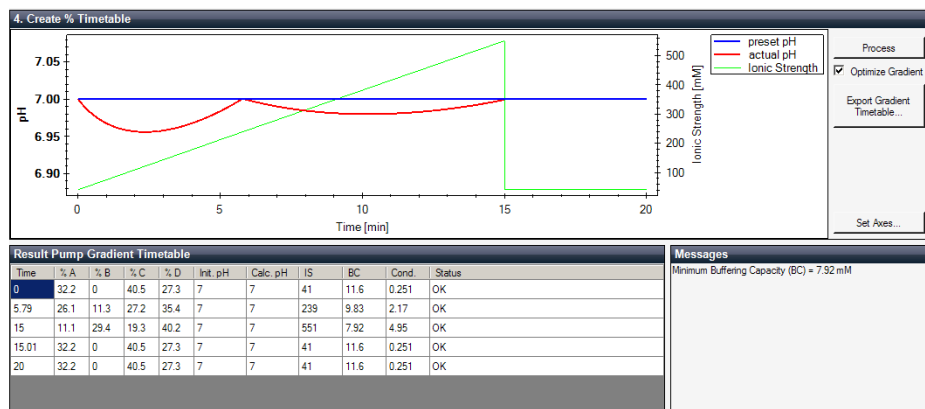
- 6 Click **Recipe** in the **Compose Stock Solutions** panel.

The **Stock Solution Recipes** dialog box is shown, containing the recipes for the stock solutions (in mg/L or mL/L). You can use these recipes to prepare the stock solutions. Bottle A contains water.



7 Click **OK** to close the **Stock Solution Recipes** dialog box.

An optimized pump gradient timetable is automatically calculated.



Short pH Gradient with Single Buffer System

pH gradients are formed by using buffers consisting of multiple buffer components. Ionic strength must be kept low in order to avoid interference with protein binding. Therefore, only bottles C and D will contain the necessary stock solutions for this experiment. In this example, your protein is stable above its pI (pH 8), therefore an anion exchange buffer is needed. The buffer system is a Tris/ HCl buffer (pKa 8.1).

- 1 In the **Select Buffer and Gradient Mode** panel, select the **Single Buffer (pH/Salt Gradient)** option.
- 2 Select the **Anion Exchange** option.
- 3 Open the dropdown menu and select **TRIS/HCl (TRIS+TRIS/HCl)** from the list of available buffers.
- 4 In the **Define Gradient Timetable** panel, enter the following gradient timetable parameters:

Time (min)	Salt	pH	Buffer
0	0	8.5	20
15	0	7.5	20
15.01	0	7.5	20
20	0	7.5	20

- 5 Click **Set** under the heading **Recommended** at the right of the **Compose Stock Solutions panel** and accept the recommended concentrations.

The data is automatically processed and the pump gradient timetable is produced.

1. Select Buffer & Gradient Mode

New... Open... Save...

Single Buffer (pH / Salt Gradient) Composite Buffer (Wide Range pH Gradient)

Cation Exchange Anion Exchange

TRIS/HCl (TRIS+TRIS/HCl)

2. Define Gradient Table

Time	Salt	pH	Buffer
0	0	8.5	20
15	0	7.5	20
15.01	0	7.5	20
20	0	7.5	20

3. Compose Stock Solutions

A: Water Recommended: 0 mM

B: NaCl [0] 0 mM

C: TRIS - HCl salt [26] 26 mM

D: TRIS [21] 21 mM

Recipe... Set

4. Create % Timetable

Process

Optimize Gradient

Export Gradient Timetable...

Set Axes...

Result Pump Gradient Timetable										
Time	% A	% B	% C	% D	Init. pH	Calc. pH	IS	BC	Cond.	Status
0	10.7	0	24.6	64.7	8.5	8.5	6.4	10.1	0.0632	OK
10.51	17.8	0	54.5	27.7	7.8	7.8	14.2	9.56	0.136	OK
15	19.9	0	63.9	16.2	7.5	7.5	16.6	6.53	0.159	OK
15.01	19.9	0	63.9	16.2	7.5	7.5	16.6	6.53	0.159	OK
20	19.9	0	63.9	16.2	7.5	7.5	16.6	6.53	0.159	OK

Messages

Minimum Buffering Capacity (BC) = 6.53 mM

You can now set up your stock solutions according to the recipe.

You can also use the single buffer mode for scouting experiments, as described in “pH Scouting with Single Buffer System” on page 38.

pH Gradient with Multi-Buffer System

pH gradients are formed by using buffers consisting of multiple buffer components. Ionic strength must be kept low in order to avoid interference with protein binding. Therefore, only bottles C and D will contain the necessary stock solutions for this experiment. The bottle containing water can be used for dilution.

Line A	Water
Line B	Not used
Line C	Acidic buffer component
Line D	Basic buffer component

- 1 In the **Select Buffer and Gradient Mode** panel, select the **Composite Buffer (Wide Range pH Gradient)** option.
- 2 Select the **Cation Exchange** option.
- 3 Open the dropdown menu and select **CEX Phosphate/Citrate, pH Range 2.8-7.1 (with salts)** from the list of available buffers.
- 4 In the **Define Gradient Timetable** panel, enter the following gradient timetable parameters:

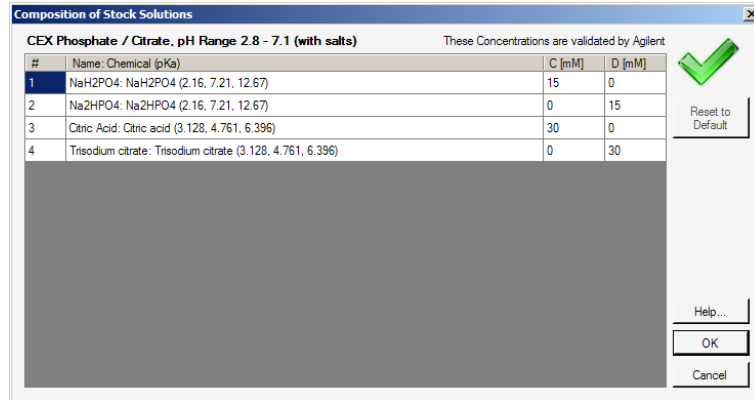
Time	pH
0	3
15	7
20	7

The screenshot shows three panels from a software interface:

- 1. Select Buffer & Gradient Mode:** Contains buttons for 'New...', 'Open...', and 'Save...'. It has radio buttons for 'Single Buffer (pH / Salt Gradient)', 'Composite Buffer (Wide Range pH Gradient)', 'Cation Exchange', and 'Anion Exchange'. A dropdown menu is set to 'CEX Phosphate / Citrate, pH Range 2.8 - 7.1'.
- 2. Define Gradient Table:** A table with columns 'Time' and 'pH'. The data rows are: (0, 3), (15, 7), (20, 7). There is a double asterisk (**) in the bottom left corner.
- 3. Compose Stock Solutions:** Lists components A, B, C, and D. Component A is 'Water' with a '%A (Dilution)' field set to 0. Component B is 'n/a'. Component C is 'pH = 2.56; IS = 18.3 mM, BC = 26.9 mM'. Component D is 'pH = 9.24; IS = 225 mM, BC = 0.267 mM'. There are 'Composition...' and 'Recipe...' buttons at the bottom.

- 5 Click **Composition** in the **Compose Stock Solutions** panel to review the composition of the stock solutions.

The **Composition of Stock Solutions** dialog box is shown, containing the suggested compositions of the components.



The green check mark indicates that these concentrations for this multi-buffer system have been validated by Agilent. However, you may change the compositions in the **Composition of Stock Solutions** dialog box and observe the change in pH range in the **Compose Stock Solutions** panel.

- 6 Click **Recipe** in the **Compose Stock Solutions** panel to display the recipes for each of the selected compositions.

The pump gradient timetable is calculated automatically and displayed.

pH Scouting with Single Buffer System

pH scouting is an approach for screening for optimum resolution of an analyte mixture using a single buffer system at different pH values.

In this example, the buffer system is a phosphate buffer (pKa 7.2) with NaCl (final concentration 500 mM). You can run a salt gradient at pH 6.5, 7 and 7.5.

First, you need to define the boundary conditions with minimum and maximum pH and minimum and maximum salt concentration to find the suitable buffer and salt stock solution that fits for the chosen pH interval.

Table 6 Salt (NaCl) gradient at pH 6.5

Time	Salt	pH	Buffer
0	0	6.5	20
15	500	6.5	20
15.01	0	6.5	20
20	0	6.5	20

Table 7 Salt (NaCl) gradient at pH 7.0

Time	Salt	pH	Buffer
0	0	7.0	20
15	500	7.0	20
15.1	0	7.0	20
20	0	7.0	20

Table 8 Salt (NaCl) gradient at pH 7.5

Time	Salt	pH	Buffer
0	0	7.5	20
15	500	7.5	20

Table 8 Salt (NaCl) gradient at pH 7.5

Time	Salt	pH	Buffer
15.01	0	7.5	20
20	0	7.5	20

1 Set up the timetable as shown in the figure:

The screenshot shows a software interface with three main panels:

- 1. Select Buffer & Gradient Mode:** Contains buttons for 'New...', 'Open...', and 'Save...'. It has radio buttons for 'Single Buffer (pH / Salt Gradient)', 'Composite Buffer (Wide Range pH Gradient)', 'Cation Exchange', and 'Anion Exchange'. A dropdown menu is set to 'Sodium Phosphate (NaH₂PO₄+Na₂HPO₄)'.
- 2. Define Gradient Table:** A table with columns 'Time', 'Salt', 'pH', and 'Buffer'. The data rows are:

Time	Salt	pH	Buffer
0	0	6.5	20
15	500	6.5	20
15.01	0	7.5	20
20	500	7.5	20
- 3. Compose Stock Solutions:** Lists components A: Water, B: NaCl, C: NaH₂PO₄, and D: Na₂HPO₄. Each has a 'Recommended' column with input fields and values: 1700, 30.5, and 37 mM respectively. A 'Recipe...' button and a 'Set' button are at the bottom.

Note that the lowest and highest salt concentrations and lowest and highest pH values are combined to define the boundary conditions for the scouting experiment.

- 2 Click **Set** in the **Recommended** section of the **Compose Stock Solutions** panel.
- 3 Note down the recommended concentrations given in the **Compose Stock Solutions** panel.

5 Ion Exchange Experiments

Generate and export a gradient time table

Generate and export a gradient time table

- 1 Click **Process** in the **Create % Timetable panel** to calculate a corrected pump gradient timetable.
- 2 Click **Export Gradient Timetable** to save the corrected gradient timetable in an appropriate file format.

The gradient time table can be saved as *.xml file format for import into the G5611A pump driver (OpenLAB CDS). From Revision A 01.04, the *.asb file format can also be used.

Import a gradient timetable into the G5611A pump driver

You can import the corrected gradient timetables into the Bio-inert pump (G5611A) driver software, which is part of OpenLAB CDS (ChemStation or EZChrom). It is supported with revision A.01.04 or later.

- 1 Go to the OpenLAB CDS ChemStation and set up and save a new method, for example, **pH6.5_salt.M**.
- 2 Go to the **Quat Pump** tab and click **Import Solvent Blending File**.
- 3 Select the file that you generated for pH 6.5 in the Buffer Advisor software (in **“Generate and export a gradient time table”** on page 40).

The timetable is filled automatically with the stored gradient values previously calculated in Buffer Advisor.

Time [min]	A [%]	B [%]	C [%]	D [%]	Flow [mL/min]	Max. Pressure Limit [bar]
0.00	37.3	0.0	49.3	13.4	0.000	600.00
0.00	37.3	0.0	49.3	13.4	---	---
4.73	29.5	9.3	40.4	20.8	---	---
15.00	11.0	29.4	31.3	28.3	---	---
15.01	43.4	0.0	14.2	42.4	---	---
15.93	38.7	5.4	10.5	45.4	---	---
17.15	31.9	12.6	8.2	47.3	---	---
20.00	15.5	29.4	5.6	49.5	---	---

- 4 Save the method.

5 Ion Exchange Experiments

Import a gradient timetable into the G5611A pump driver

5 Generate new methods for pH 7.0 and pH 7.5.

You have now generated three methods in OpenLAB CDS ChemStation with the respective gradient timetables. **Figure 8** on page 42 shows an example chromatogram of phosphate buffer salt gradients run at different pH.

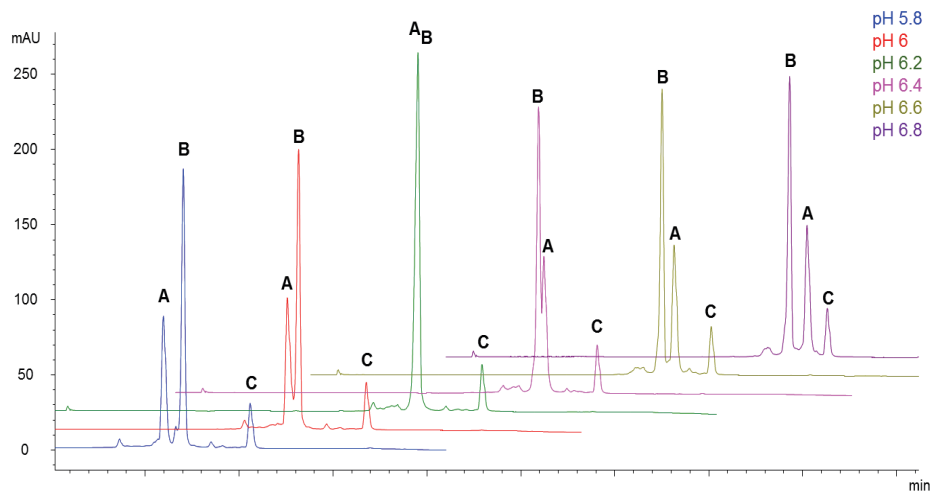


Figure 8 pH scouting experiment using cytochrome C, lysozyme and RNase.



6 Troubleshooting and Tips

Messages shown by buffer advisor 44

Essential measurement practices 46

This chapter explains the messages that the Agilent Buffer Advisor provides, and gives you some important information on how to avoid problems.



Messages shown by buffer advisor

Messages in Buffer Advisor software give you hints on how to optimize parameters in the software, such as stock solution concentration or gradient timetable entries for best separation conditions.

Perhaps the buffering capacity (BC) of the buffer is very low, for example, less than 3 mM, and the buffer might not buffer any more, and therefore the entered pH cannot be maintained.

Another example is that a liquid chromatograph's proportioning valve cannot mix one of the solutions precisely if the added percentage of this solution is too low. For best performance, it is recommended always to mix more than 5% per channel.

Here are some parameters that can be modified:

- *Buffer (pH range)*: Select a different buffer (system) closer to the pH value you entered in the gradient timetable.
- *Salt concentration (around 500 - 1000 mM)*: Ionic strength can influence pH of buffers by lowering buffering capacity. Lower the salt concentration in case of errors and see if the buffering capacity increases.
- *pH*: When you want to keep the same buffer you already selected, change your desired pH in the gradient timetable to ± 0.5 of pKa of buffer.
- *Buffer concentration (around 10 - 50 mM)*: Increase buffer concentration when your buffering capacity is low or when you get a message that pump mixing is less than 5%.
- *Stock solutions*: Decrease concentration of stock solutions when the 5% warning is shown.

Table 9 List of messages and proposed actions

Message	Action
Calculation process detected database error.	Stay within pH range of selected buffer, lower salt concentration or increase buffer concentration.
Unexpected calculation error ("{0}").	Stay within pH range of selected buffer, lower salt concentration or increase buffer concentration.

Table 9 List of messages and proposed actions

Message	Action
Timetable input cannot be calculated with selected buffer. Select different buffer or change timetable.	Select different buffer or change timetable entries to adjust gradient. Adjust parameters as shown above.
Minimum Buffering Capacity (BC) = {0} mM	Only take action when other messages are shown and try to keep it above 3 mM.
Increase mixing per channel > 5% by editing gradient timetable. Choose pH near PKa of buffer, increase buffer concentration or lower salt concentration.	Stock solutions are highly concentrated, leading to a low mixing proportion (less than 5%) of the quaternary pump. Lower stock solutions concentration or increase buffer concentration.
pH difference larger than set threshold {0} detected in "pH Difference" plot.	The pH entered in the gradient table is too far away from the buffer's pKa. Choose pH near pKa, or increase buffer concentration.
Maximum number of lines in result pump gradient timetable reached (maximum is {0} lines). Check pH differences if acceptable.	The pH entered in the gradient table might be too far away from the buffer's pKa so that the corrected gradient timetable contains a lot of mixing steps. Try to stay close to the pKa of the buffer.
Calculation of one or more result pump gradient timetable lines ended with error. Check their status.	Check status line in pump timetable, then adjust parameters.
Stock solution concentration for '{0}' in Bottle {1} exceeded concentration of stock liquid. Maximum concentration is {2} mM.	Final buffer concentration is higher than stock solution concentration. Decrease buffer concentration or increase stock solution concentration.
'{0}' in Bottle {1} may be insoluble in given stock solution concentration. Maximum solubility is expected around {2} mM.	Consider different stock solution concentrations.

Essential measurement practices

Storage and Handling

- Handle and store all reagents according to MSDS and the instructions on the label of the individual box.
- Keep all reagent and reagent mixes refrigerated at 4 °C when not in use.
- Allow all reagents and samples to equilibrate to room temperature for 30 minutes before use.
- Use clean bottles and material for buffers and use fresh solvents to avoid microbial growth.

Preparation of Stock Solution Buffers and Salt Solutions

- The buffering substance should have the same charge as the exchanger.
- For the adjustment of pH, the same counter-ion should be used as for elution (e.g. HCl/NaCl)
- Buffer contaminants may produce “ghost” peaks. Use highly purified buffer salts.
- Always use a buffering agent that has a pKa within ± 0.5 pH units of the operating pH. Otherwise, there is a risk of dramatic pH fluctuation due to the limited buffering capacity of the buffering agent.
- pH varies with the temperature. Tris buffers are particularly temperature sensitive. Adjust pH value after buffers have been equilibrated to the desired running temperature.
- Keep in mind that especially strong bases, for example, NaOH, take CO₂ from air to build carbonate, leading to a change of pH over time.
- Temperature can influence pH. All calibrations and measurements should be performed at the same temperature.
- Note that pH electrodes age and should be replaced on regular basis (i.e. once per year) and/or when pH slope and zero voltage offset are out of recommended range specified by manufacturer.
- Always calibrate your electrode before measurement of pH within the appropriate pH of calibration solutions. Do not use expired calibration solutions and do not reuse them.

- If measuring pH at high ionic strength make sure that your pH electrode is suited for this task.

Mixing Properties of Agilent 1260 Infinity LC Quaternary Pump

We have taken great care that the pump properties of the 1260 Infinity Bio-inert LC system are taken into account for calculation of optimum conditions of the gradients.

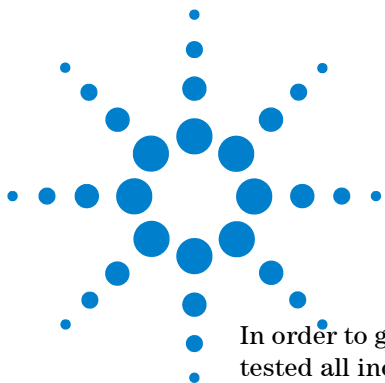
For example, the channels defined for separation must always be A for Water, B for salt and C and D for acidic and basic component.

Be sure to flush the channels properly after use. High concentrations of salt can result in the formation of crystals that can impede the function of the proportioning valve.

- For highest pH precision:
 - Set primary channel in pump driver interface to “A” (accessible via the advanced tab).
 - Make sure that the percentage of acidic and basic solution is always above 5% of the total mobile phase in order to ensure highest gradient precision.

6 Troubleshooting and Tips

Essential measurement practices



7

Validation Tests and Specifications

In order to guarantee high accuracy of the resulting pH of the mobile phase we tested all indicated buffers under the following conditions:

Instrumentation

Bio-inert quaternary pump (G5611A) connected to a thermostatted column compartment (G1316C with bio-inert heat exchanger G5616-60050), Diode-array detector (G1315C with bio-inert 10 mm path length flow cell G5615-60022) and bio-inert analytical fraction collector (G5664A).

Method for salt gradient mode buffers

Each buffer has been tested at

- three different buffer concentrations (10, 50, 100 mM)
- three different pH values (approximately ± 0.5 pH units distance to published pKa value)
- at least three (up to five) different salt concentrations ranging from 0 to 1000 mM NaCl (5 min salt step gradient)

1 mL fractions were taken over the whole run and pH values were measured offline using an appropriately calibrated pH sensor (Mettler Toledo Inlab micro, Schwerzenbach, Switzerland) connected to a pH meter (Schott handylab pH/LF 12, Mainz, Germany).

All tests were carried out at 23 °C. If the separation in your application takes place at a different temperature, deviations from the set pH can occur

Method for pH gradient multicomponent buffers

After setting up the two bottles with the indicated constituents and concentrations, a pH step gradient method was performed. Each pH step lasted for 5 min, pH difference between the steps was 0.5 – 1.0 pH units.



7 Validation Tests and Specifications

Essential measurement practices

1 mL fractions were taken over the whole run and pH values were determined according to the “[Method for salt gradient mode buffers](#)” on page 49.

Stock solutions

Buffer Advisor was used to simulate validation runs in order to identify the minimum number of buffer compound stock solutions necessary to run all validation runs for the corresponding buffer (typically 2 to 3 concentrations of acidic and basic buffer solutions, respectively). Care was taken to maintain acidic and basic channels (C and D) above 5% for highest mixing precision (see “[Essential measurement practices](#)” on page 46).

The maximum deviation from set pH in all buffers tested is +/- 0.2 pH units.¹

Validated salt gradient buffer list

The following buffers have been tested according to the procedure stated above.

Cation exchange buffers		Anion exchange buffers	
Buffer Name	pH tested	Buffer Name	pH tested
Na-phosphate	6.2, 7.0, 7.7	Piperazine	5.2, 5.7, 6.2, 9.3, 9.8, 10.3
Na-citrate	4.3, 4.8, 5.3, 5.9, 6.4, 6.9	Bis-Tris	6.0, 6.5, 7.0
Formic acid	3.2, 3.7, 4.2	Triethanolamine	7.3, 7.8, 8.3
Lactic acid	3.4, 3.9, 4.4	Tris	7.6, 8.2, 8.6
Acetic acid	4.3, 4.8, 5.3	Diethanolamine	8.4, 8.9, 9.4
Succinic acid		Ammonia	8.6, 9.1, 9.6
Malonic acid	5.2, 5.7, 6.2	Ethanolamine	9.0, 9.5, 10.0
MES	5.6, 6.1, 6.6	Piperidine	
Maleic acid	5.8, 6.3, 6.8		
MOPS	6.7, 7.2, 7.7		
HEPES	7.1, 7.6, 8.1		

¹ Some buffers or combination of buffers were not tested and therefore may deviate from the above specifications.

Cation exchange buffers		Anion exchange buffers	
Buffer Name	pH tested	Buffer Name	pH tested
TAPS	8.1, 8.6, 9.1		
Na-bicarbonate	9.3, 9.8, 10.2		

Validated multi-component buffer list

Multi-component buffer CEX Phosphate/Citrate, pH Range 3 – 7 (with salts) was tested at following concentrations and conditions:

	Bottle C	Bottle D
Na ₂ HPO ₄		15 mM
NaH ₂ PO ₄	15mM	
citric acid	30 mM	
trisodium citrate		30 mM
pH calculated	2.55	9.25
pH measured	2.50	9.04

A pH gradient was run from pH 3.0 to pH 7.5.

Multi-component buffer AEX, pH Range 6 – 11 (with salts) was tested at following concentrations and conditions:

	Bottle C	Bottle D
Bis-Tris		60 mM
Bis-Tris-HCl salt	60 mM	
Tris		30 mM
Tris-HCl salt	30 mM	
Ethanolamine		40 mM
Ethanolamine-HCl salt	40 mM	
Piperidine		35 mM

7 Validation Tests and Specifications

Essential measurement practices

	Bottle C	Bottle D
Piperidine-HCl salt	35 mM	
pH calculated	3.94	11.79
pH measured	3.97	11.80

A pH gradient was run from pH 6.0 to pH 11.0.

Index

A

anion exchange chromatography 18
anion exchange 10, 20

B

buffer concentration 6, 44
buffer selection 26
buffer system 26
buffer 11, 15, 18, 20, 44
buffering capacity 44

C

cation exchange chromatography 18
cation exchange 10, 20
check mark 28
chromatofocusing 14
composite buffer mode 25
composite buffer 26, 36
counter-ion 18

D

demo mode 24

G

gradient timetable 26

I

i strength 6
installation 24
ion exchange chromatography 6, 10
ionic strength 11, 20, 44
isoelectric point 10

L

license key 24

P

pH gradient 6, 10, 14, 25, 34, 36
pH scouting 6, 38
pH 10, 20, 44
pI 10
pKa 10, 20, 27
pump timetable 6

R

recipe 26, 28

S

salt concentration 6, 44
salt gradient 6, 10, 12, 25, 32
single buffer mode 25
single buffer 26, 32, 34, 38
stock solution 7, 26, 44

T

timetable 6

www.agilent.com

In This Book

This manual describes the Agilent Buffer Advisor software and its use. It gives background information on ion exchange chromatography, and describes typical workflows for defining stock buffer solutions for ion exchange experiments.

© Agilent Technologies 2012

Printed in Germany
5/12



G5617-90000



Agilent Technologies