

CX-1 pH Gradient Buffer
QuickStart

- 083273 CX-1 pH Gradient Buffer A (pH 5.6), 125 mL
- 085346 CX-1 pH Gradient Buffer A (pH 5.6), 250 mL
- 083275 CX-1 pH Gradient Buffer B (pH 10.2), 125 mL
- 085348 CX-1 pH Gradient Buffer B (pH 10.2), 250 mL

Lot Number YYMMDD

Expiration of Certification
Month YYYY

1. Overview

- CX-1 pH Gradient Buffers are designed to generate a linear pH gradient with cation exchange columns for the characterization of MAb charge variants and other proteins.
- This QuickStart guide will provide step-by-step instructions to prepare the pH Gradient Buffer concentrates.
- For more detailed information regarding the use of pH gradient buffers, please consult the pH Gradient Buffer User Manual, Document No. 065534, at <http://www.thermoscientific.com/dionex>.

2. Main Features

- Generates robust and reproducible pH gradients.
- Ready to use with existing LC columns and systems without the need for time consuming mobile phase adjustments.

3. Directions for Use

- Dilute the 10X pH Gradient Buffer A or B 10-fold using either weight/weight or volume/volume with Type 1 reagent grade deionized water with a specific resistance of 18.2 Megohm-cm or greater filtered through a 0.2 µm filter.
- Set the UV wavelength to 280 nm or above.
- To generate a pH gradient from 5.6 to 10.2 set the gradient program from 100% A to 100% B over 30 minutes for MAbPac SCX-10, 10 µm, 4 × 250 mm column. For other columns, adjust the gradient time based on the column dimension.

NOTES:

- Store the 10X pH Gradient Buffer A and Buffer B between 2-8 °C.
- If the buffers are frozen, equilibrate to room temperature and mix well prior to diluting.
- DO NOT dilute Buffer A or Buffer B more than 10-fold. Over dilution will reduce the overall buffering capacity.
- To achieve optimal chromatographic separation, it is recommended to use Thermo Fisher Scientific ProPac WCX-10 and MAbPac SCX-10 columns.

CERTIFICATE OF ANALYSIS

The CX-1 pH Gradient Buffers were developed, formulated, and tested for use with the ProPac WCX-10 and MAbPac SCX-10 columns. The preparation and analyses, for pH, conductivity and UV absorption, of this lot of pH Gradient Buffers was performed with extreme care by Thermo Scientific Consumables Manufacturing Department in Sunnyvale California to ensure this lot meets the specifications stated below.

| Parameter | CX-1 pH Gradient Buffer A | CX-1 pH Gradient Buffer B |
|-------------------------------|---------------------------|---------------------------|
| pH | 5.6 ± 0.2 | 10.2 ± 0.2 |
| Conductivity (mS/cm) | 26.0 ± 2.0 | 40.0 ± 2.0 |
| UV absorption @ 280 nm | ≤ 0.01 | ≤ 0.01 |

A full Certificate of Analysis can be made available upon request by contacting techsupport@dionex.com

Thermo Scientific offers a 3-month warranty on this product. Tests have been completed to ensure that concentration remains within the certified values throughout the product shelf life.