

Profinia™ Bottled Buffers

Catalog No.

620-0205, 620-0206, 620-0207, 620-0208, 620-0209, 620-0210, 620-0211, 620-0212, 620-0213, 620-0214, 620-0215, 620-0216, 620-0217, 620-0218, 620-0219

Description

The Profinia protein purification system provides buffers and reagents for the purification of affinity-tagged proteins using native IMAC, denaturing IMAC, GST, or desalting methods. Bottles of buffers and reagents are available individually for the convenience of customizing buffer packs for use with the custom methods on the Profinia instrument.

Storage

Store between 4 and 22°C

Instructions for Use

All Profinia buffers are supplied as concentrates and the concentration of each solution should be programmed into the custom methods of the instrument. The denaturing IMAC buffers and GST elution buffer require the addition of urea (620-0200 and 620-0201) or glutathione packs (620-0202) prior to use. Unused solutions can be stored at –20°C for up to 6 months and thawed once prior to use.

Shelf Life

The buffers are stable for at least one year from date of receipt. Refer to the specific expiration date on the product label.

Formulations

| Item # | Description | Supplied As | Formulations (of concentrates) |
|----------|--------------------------------|-------------|--|
| 620-0205 | Native IMAC Lysis Buffer | 2x | 600 mM KCl, 100 mM KH ₂ PO ₄ , 10 mM Imidazole, pH 8.0 |
| 620-0206 | Native IMAC Wash Buffer-1 | 2x | 600 mM KCl, 100 mM KH ₂ PO ₄ , 10 mM Imidazole, pH 8.0 |
| 620-0207 | Native IMAC Wash Buffer-2 | 2x | 600 mM KCl, 100 mM KH ₂ PO ₄ , 20 mM Imidazole, pH 8.0 |
| 620-0208 | Native IMAC Elution Buffer | 2x | 600 mM KCl, 100 mM KH ₂ PO ₄ , 500 mM Imidazole, pH 8.0 |
| 620-0209 | Denaturing IMAC Lysis Buffer | 1.4x | 420 mM KCl, 70 mM KH ₂ PO ₄ , 7 mM Imidazole, pH 8.0 |
| 620-0210 | Denaturing IMAC Wash Buffer-1 | 1.4x | 420 mM KCl, 70 mM KH ₂ PO ₄ , 7 mM Imidazole, pH 8.0 |
| 620-0211 | Denaturing IMAC Wash Buffer-2 | 1.4x | 420 mM KCl, 70 mM KH ₂ PO ₄ , 14 mM Imidazole, pH 8.0 |
| 620-0212 | Denaturing IMAC Elution Buffer | 1.4x | 420 mM KCl, 70 mM KH ₂ PO ₄ , 350 mM Imidazole, pH 8.0 |
| 620-0213 | GST Lysis Buffer | 2x | 300 mM NaCl, 20 mM Na ₂ HPO ₄ , 10 mM EDTA, pH 7.4 |
| 620-0214 | GST Wash Buffer | 2x | 300 mM NaCl, 20 mM Na ₂ HPO ₄ , 10 mM EDTA, pH 7.4 |
| 620-0215 | GST Elution Buffer | 2x | 40 mM Glutathione, 200 mM Tris, 10 mM EDTA, pH 8.0 |
| 620-0216 | Desalting Buffer | 5x | 685 mM NaCl, 13.5 mM KCl, 21.5 mM Na ₂ HPO ₄ , 40.5 mM KH ₂ PO ₄ , pH 7.0 (pH 7.4 upon dilution) |
| 620-0217 | Cleaning Solution 1 | 2x | 1,000 mM NaCl, 100 mM Tris, pH 8.0 |
| 620-0218 | Cleaning Solution 2 | 4x | 2,000 mM NaCl, 400 mM NaOAc, pH 4.5 |
| 620-0219 | Storage Solution | 2x | 4% C ₆ H ₅ CH ₂ OH (Benzyl Alcohol) |