
**Profinia™ Denaturing
IMAC Urea Pack**

45 g, 90 g

**Catalog Numbers
620-0200, 620-0201**

BIO-RAD

Specifications

Quantity	45 g, 90 g
Description	Profinia urea packs contain high-purity chromatography grade urea which is ideal for solubilizing and denaturing proteins in <i>E. coli</i> extracts.
Storage	Store between 4°C and 22°C
Instructions for use	Add the entire contents of the 90 g pack to the 176 ml of Profinia Denaturing IMAC Lysis Buffer (#620-0209) or Profinia Denaturing IMAC Wash Buffer 1 (#620-0210) to make 250 ml of a 6 M urea solution. Add the entire contents of the 45 g pack to the 88 ml of Profinia Denaturing IMAC Wash Buffer 2 (#620-0211) or Profinia Denaturing IMAC

Elution Buffer (#620-0212) to make 125 ml of a 6 M urea solution. Stir to dissolve, and filter through a 0.45 μm filter to remove any particulates. Fresh urea solutions should be prepared and used the day of chromatography. Unused solutions can be stored at 4°C and be used within 7 days, or stored at -20°C for up to 6 months and thawed once prior to use.

Shelf life

The urea packs are stable for at least 2 years from date of receipt. Refer to the specific expiration date on the product label.

