
Glutathione Pack

1.23 g

**Catalog Number
620-0202**

BIO-RAD

Specifications

Quantity	1.23 g
Description	The glutathione pack contains high-purity reduced glutathione, which is ideal for eluting glutathione S-transferase (GST) fusion proteins from immobilized glutathione resins.
Storage	Store at 4°C
Instructions for use	Add the entire contents of the 1.23 g pack to the 100 ml of GST Elution Buffer (catalog #620-0215) to generate a 40 mM glutathione 2x solution (final working concentration of glutathione is 20 mM). Stir to dissolve and filter through a 0.45 µm filter to remove any particulates. Fresh glutathione

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 glutathione packs are stable for at least 1 year from date of receipt. Refer to the specific expiration date on the product label.

