

Data Sheet

PARP9, GST-Tag, His-Tag

Human, recombinant, N-terminal GST-tag,
C-terminal His-tag

Catalog #: 80509

Lot#: 200716-A

Conc.: 0.32 mg/ml

Description: Human PARP9, GenBank Accession No. NM_031458, a.a. 1-854(end) with N-terminal GST-tag and C-terminal His-tag, expressed in a Baculovirus infected Sf9 cell expression system. MW = 123 kDa.

Formulated in: 40 mM Tris-HCl, 110 mM NaCl, 2.2 mM KCl, 3 mM DTT, 20% Glycerol and 20 mM Glutathione.

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles.

References:

1. Hakmé A., *et al. Dev. Dyn.* 2008. Jan; **237(1)**:209-15.
2. Moyle PM, Muir TW. *J Am Chem Soc.* 2010. Nov 17; **132(45)**:15878-80.

Quality Assurance

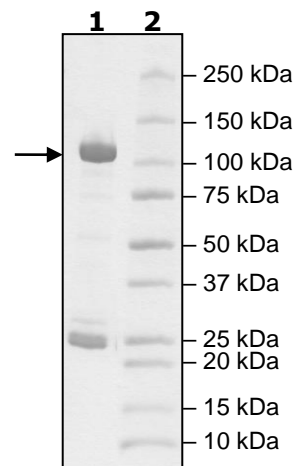
4-20% SDS-PAGE Coomassie staining

Lane 1:
3 μg PARP9

Lane 2:
Protein Marker

MW: 123 kDa

Purity: 58%



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com