

Data Sheet

PARP1, Flag-Avi-tag

Human, recombinant, N-terminal FLAG-Avi-tag

Catalog #: 80521

Lot #: 150715-1 **Conc.:** 0.70 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, 0.2mM TCEP, 80 µg/ml FLAG peptide

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

References:

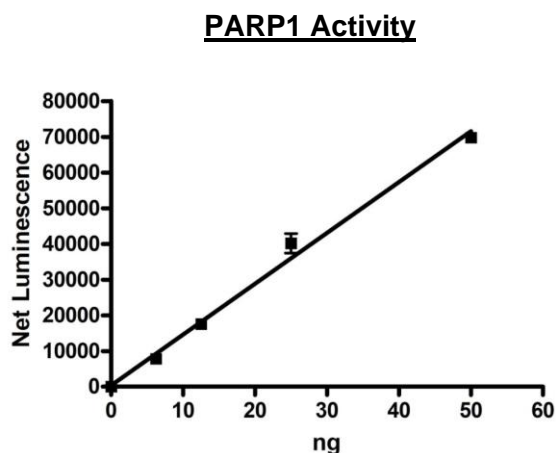
1. Aguilar-Quesada R, *et al.*, *Curr Med Chem.* 2007;**14(11)**:1179-87.
2. Pacher P. and Szabó C. *Cardiovasc Drug Rev.* 2007 Fall;**25(3)**:235-60.

Description: Human PARP1 (GenBank Accession No. NM_001618), full length with N-terminal FLAG-Avi-tag™, MW = 117 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

Assay Conditions: Enzyme reaction is conducted using PARP Assay Kit (BPS Bioscience, #80551) at room temperature for 1 hour.

Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance



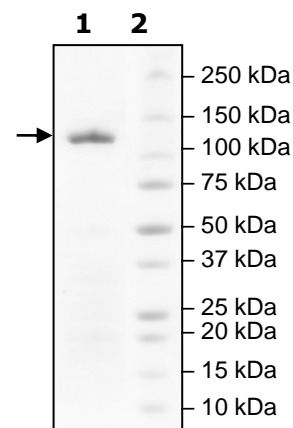
4-20% SDS-PAGE Coomassie staining

Lane 1:
3.8 µg PARP1

Lane 2:
Protein Marker

Purity: > 90%

MW: 117 kDa



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