

## Data Sheet

### PARP1, GST-tag

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

**Catalog #:** 80521

**Lot #:** 150715-1

**Conc.:** 0.70

**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. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).*

**References:**

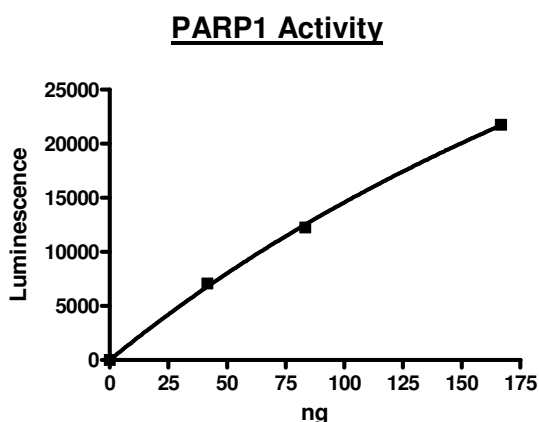
1. Aguilar-Quesada R, et al., *Curr Med Chem.* 2007;**14(11)**:1179-87.
2. Pacher P. and Szabó C. *Cardiovasc Drug Rev.* 2007 Fall;**25(3)**:235-60.

**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 BPS PARP Assay Kit (Catalog Number 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

