

Data Sheet

PARP10, FLAG-Strep-Tag

Human, Recombinant, N-terminal FLAG-Tag,
 C-terminal Strep-Tag

Catalog #: 80522

Lot #: 161111 **Conc.:** 0.56 mg/ml

Description: Human PARP10, GenBank Accession No. NM_032789, with N-terminal FLAG-Tag and C-terminal Strep-tag, a.a. 805-1025, MW = 27 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween20, 20% glycerol, 80 µg/ml flag peptide

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

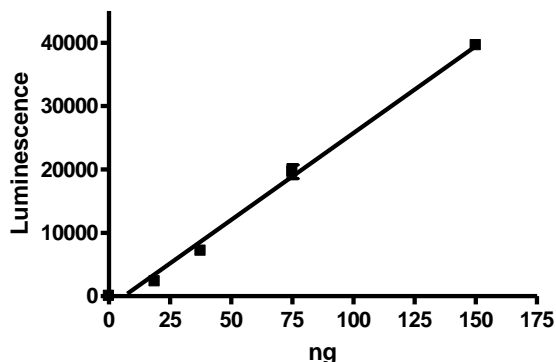
Stability: At least 6 months at -80°C. Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

References:

1. Kleine, H., et al., *Molecular Cell*. 2007; **32(1)**:57-69.
2. Chou, H., et al., *Journal of Biological Chemistry*, 2006 Feb; **281**:15201-15207.

Quality Assurance

Specific Activity



4-20% SDS-PAGE Coomassie staining

Lane 1:
 PARP10
 1 µg loaded

Lane 2:
 Protein Marker

MW: 27 kDa
Purity: ≥86%

