

## Data Sheet

### **TNKS2(667-end) (PARP5B), GST-tag**

Human, recombinant, N-terminal GST-tag  
**Catalog #:** 80505  
**Lot #:** 130314-G2      **Conc.:** 0.28 mg/ml

**Formulated in:** 40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween, 20% Glycerol, and 3 mM DTT.

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

**References:**

1. Cook, B.D. *et al.* (2002). *Mol. Cell. Biol.* **22**(1): 332–42.
2. Kaminker PG, *et al.* (2001). *J. Biol. Chem.* **276** (38): 35891–9.
3. Sidorova N, *et al.* (2006). *Cell Tissue Res.* **323** (1): 137–45.

**Description:**

TNKS2 (TRF1-interacting ankyrin-related ADP-ribose polymerase 2), also known as PARP5B, TNKL, or tankyrase-2 (GenBank Accession No. NM\_NM\_025235), a.a. 667-1166(end), with N-terminal GST-tag, MW=80.5 kDa, expressed in Sf9 cells via a baculovirus expression system.

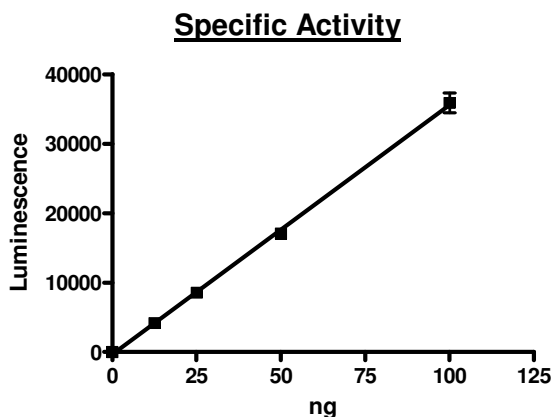
**Assay Conditions:**

Assay was performed following the BPS Bioscience TNKS2 Histone Ribosylation Assay Kit (Biotin-labeled NAD+) protocol, Catalog # 80578.

**Application:**

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

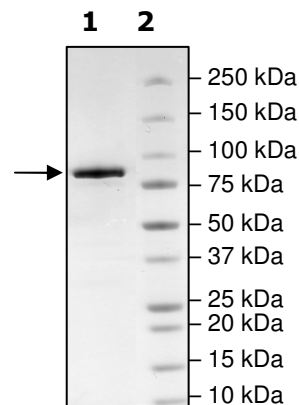
## **Quality Assurance**



**4-20% SDS-PAGE  
Coomassie staining**

**Lane 1:**  
2.5  $\mu\text{g}$  TNKS2  
**Lane 2:**  
Protein Marker

**MW:** 80.5 kDa  
**Purity:**  $\geq 97\%$



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

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**  
 Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)  
 Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)