

## Data Sheet

### DPP7, His-Tag

Human, Recombinant, C-terminal His tag

**Catalog #:** 80070

**Lot#:** 111207-2      **Conc.:** 0.07 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 200 mM imidazole, and 20% glycerol.

**Stability:** >6 months at -80°C

**References:**

1. Shreder K.R. et al., Bioorg Med Chem Lett. 2005 Oct 1;15(19):4256-60.
2. Maes M.B. et al., Biochem J. 2005 Mar 1;386(Pt 2):315-24.

**Description:**

Human DPP7 (GenBank Accession No. NM\_013379), (amino acids 26-end) with C-terminal His tag, MW=53 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

**Specific Activity:** 1880 pmol/min/μg

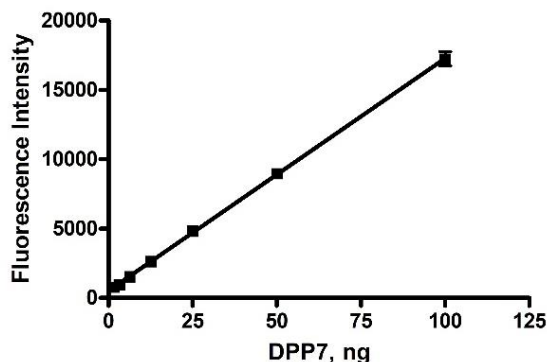
**Assay Condition:** Test wells contained 0.01M Tris-HCl pH7.4, 0.01M MgCl<sub>2</sub>, 0.05% Tween-20, 5 μM DPP substrate 1, and various amounts of DPP7(0-200ng). Standard wells contained 0.01M Tris-HCl pH7.4, 0.01M MgCl<sub>2</sub>, 0.05% Tween-20, and various amounts of Fluorescent AMC standard (0.01 μM-1.25 μM). Incubated for 10 minutes at room temperature. Fluorescence intensity was measure at excitation 360 nm and emission 460 nm.

**Application:**

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

### Quality Assurance

**Specific Activity**

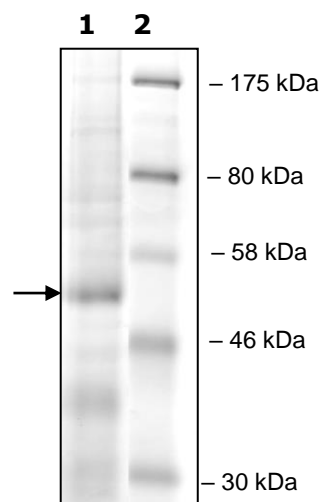


**10% SDS-PAGE  
Coomassie staining**

**Lane 1:**  
1.4 μg DPP7

**Lane 2:**  
Protein Marker  
BioLabs (#P7708L)

**MW:** 53 kDa.  
**Purity:** ≥47%



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)