

## Data Sheet

### HDAC5, GST-Tag, His-Tag

Human, Recombinant, N-terminal GST-tag,  
 C-terminal His-tag

**Catalog #:** 100102

**Lot #:** 180709-2      **Conc.:** 0.32 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 16mM Glutathione, 20% glycerol.

**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:**

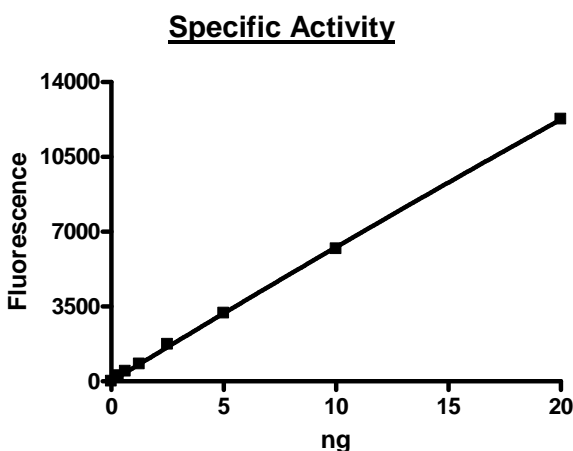
1. Spiegelberg, B.D. and Hamm, H.E. (2005) *J. Biol. Chem.* **280**, 41769-41776
2. Waltregny et al., (2004) *J. Neural Transm.* **48**, 273-290.

**Description:** Human HDAC5, also known as histone deacetylase 5, GenBank Accession No. NM\_005474, a.a. 1-1122 (end) with N-terminal GST-tag and C-terminal His-tag, expressed in a HEK293 expression system. MW=149 kDa.

**Specific Activity:**  $\geq 1000$  pmol/min/ $\mu\text{g}$ .  
 Assay Conditions: 25 mM Tris/HCl, pH 8.0, 137 mM NaCl, 2.7 mM KCl, 1 mM  $\text{MgCl}_2$ , and 0.1 mg/ml BSA, 20  $\mu\text{M}$  BPS HDAC substrate 1 (Catalog #50040), and HDAC5 (0.3125 – 20 ng).  
 Incubation condition: 30 min at  $37^{\circ}\text{C}$  followed by developing for 15 min at room temperature.

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

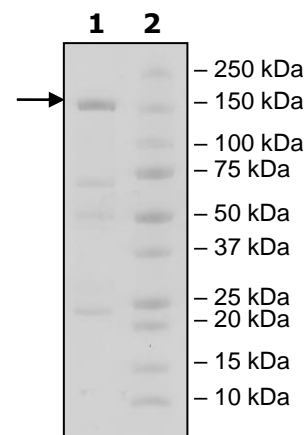
## Quality Assurance



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

**Lane 1:**  
 2  $\mu\text{g}$  HDAC5  
**Lane 2:**  
 Protein Marker

**MW:** 149 kDa  
**Purity:**  $\geq 55\%$



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)