

## Data Sheet

### HDAC 5 Full Length

Human, recombinant, N-terminal GST tag  
Catalog #: 50045  
Lot #: 130620      Conc.: 0.04 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 18 mM Glutathione and 20% Glycerol.

**Stability:** >6 months at -80 °C

#### 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.

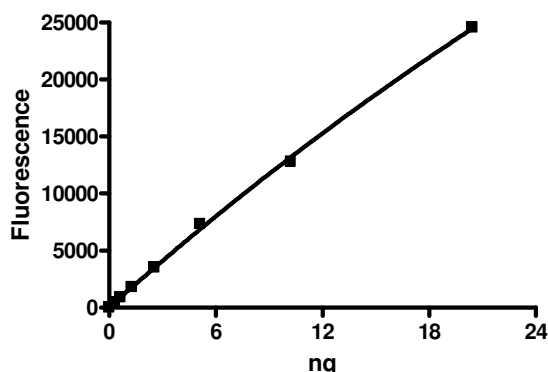
**Description:** Human HDAC5 (GenBank Accession No. NM\_005474), full length with N-terminal GST tag, MW= 150 kDa, expressed in baculovirus expression system.

**Specific Activity:** 940 pmol/min/μg.  
Assay condition: 25 mM Tris/HCl, pH 8.0, 137 mM NaCl, 2.7 mM KCl, 1 mM MgCl<sub>2</sub>, and 0.1 mg/ml BSA, 20 μM BPS HDAC substrate 1 (Catalog #50040), and HDAC5 (0.8-50 ng). Incubation condition: 30 min at 37°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

### Specific Activity

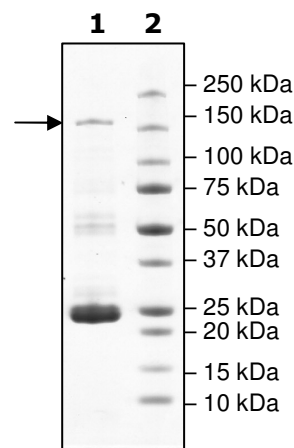


### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
1 μg HDAC5

**Lane 2:**  
Protein Marker

**MW:** 150 kDa  
**Purity:** ≥8%



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)