



11526 Sorrento Valley Rd SteA2  
San Diego Ca 92121  
Tel: 1.858.202.1401  
1.858.829.3082  
Fax: 1.858.259.5734  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## Data Sheet

### **HDAC 1**

human, recombinant, C-terminal GST tag  
Catalog #: 50001

**Formulated in:** 25 mM Tris-HCl, pH 8.0, 130 mM NaCl, 0.05% Tween-20, 20 mM glutathione and 10% glycerol.

**Stability:** >6 months at -80°C

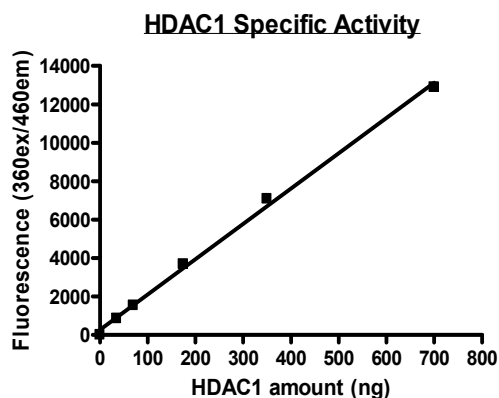
**References:**

1. Rikimaru, T. et al., *Oncology* 72 (1-2), 69-74 (2007).
2. Ishihama, K. et al., *J Clin Pathol* 60(11), 1205-10 (2007).

**Description:** Human HDAC1 (GenBank Accession No. NM\_004964), full length with C-terminal GST tag, MW= 79.9 kDa, expressed in baculovirus expression system.

**Specific Activity:** 2000 units. 1 unit is defined as the amount of enzyme that will deacetylate 1 pmol HDAC peptide substrate per minute at pH 7.4 and 37°C. Assay condition : 25 mM Tris/Cl, pH8.0, 137 mM NaCl, 2.7 mM KCl, 1 mM MgCl<sub>2</sub>, and 0.1 mg/ml BSA, 20 µM BPS HDAC substrate (Catalog number 50037), and 4 ng/µl HDAC1. Incubation condition: 30 min at 37°C.

### **Quality Assurance**

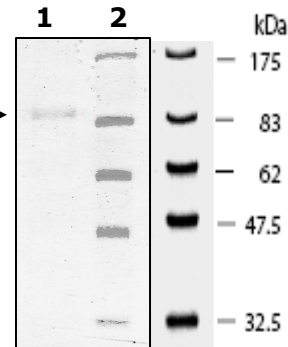


**10% SDS-PAGE**  
**Coomassie staining**

**Lane 1:**  
1 µg HDAC1 →

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

**MW:** 79.9 kDa.



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

To place your order, please contact us by Phone **1.858.202.1401** or **1.858.829.3082**, Fax **1.858.259.5734**  
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)



11526 Sorrento Valley Rd SteA2  
San Diego Ca 92121  
Tel: 1.858.202.1401  
1.858.829.3082  
Fax: 1.858.259.5734  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## Assay Protocol

**Material:** Assay buffer ( BPS catalog number 50031); Assay developer (BPS catalog number 50030); HDAC Substrate (BPS number 50037)

**Step 1:** adding all reaction mixture to a low binding NUNC black plate (VWR catalog number 62408-936)

35  $\mu$ l of HDAC assay buffer (BPS catalog number 50031)  
5  $\mu$ l of 1 mg/ml BSA  
5  $\mu$ l of 200 uM substrate (BPS catalog number 50037)  
5.0  $\mu$ l of HDAC1 (2 U/ $\mu$ l)

Always add HDAC1 at the last.

Incubate at 37 °C for 30 min.

**Setp2:** stop the reaction

add 50  $\mu$ l of HDAC assay developer (2x) (BPS catalog number 50030) and incubate the plate at room temperature for 15 min.

**Step 3:** read sample in a microtiter-plate reading fluorimeter capable of excitation at a wavelength in the range 350-380 nm and detection of emitted light in the range 440-460 nm.

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

To place your order, please contact us by Phone **1.858.202.1401** or **1.858.829.3082**, Fax **1.858.259.5734**

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)