



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 2**

Human, recombinant, C-terminal His tag  
Catalog #: 50002

**Formulated in:** 25 mM Tris-HCl, pH 8.0, 138 mM NaCl, 0.05% Tween-20 and 10% glycerol.

**Stability:** >6 months at -80°C

#### **References:**

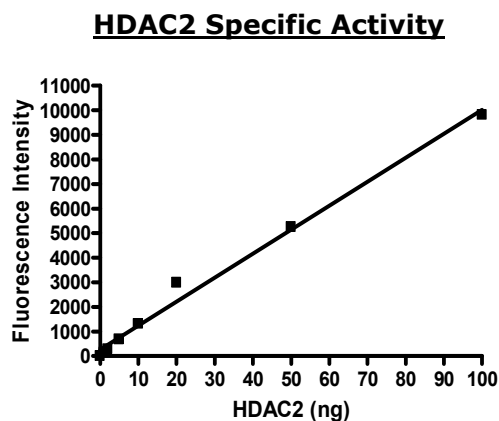
1. Yang, W.M. et al., Proc. Natl. Acad. Sci. U.S.A. 93 (23), 12845-12850 (1996)
2. Mungall, A.J. et al., Nature 425 (6960), 805-811 (2003)

**Description:** Human HDAC2 (GenBank Accession No. Q92769), full length with C-terminal His tag, MW= 60 kDa, expressed in baculovirus expression system.

**Specific Activity:** 450 pmol/min/μg. Assay condition: 50 mM of Tris-HCl (pH 8.0), 137 mM of NaCl, 2.7 mM of KCl, 1 mM of MgCl<sub>2</sub>, 5 μg BSA, 20 μM substrate (BPS catalog #: 50037), 37°C, 30 min.

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

### **Quality Assurance**

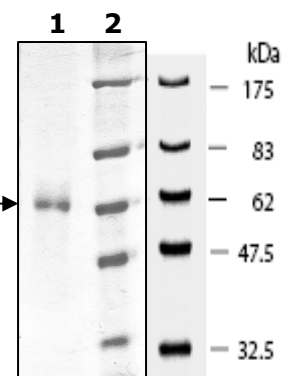


#### **10% SDS-PAGE Coomassie staining**

**Lane 1:**  
2 μg HDAC2

**Lane 2:**  
Protein Marker

**MW:** 60 kDa.  
**Purity:** >95%



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 50032)

### **Step 1: adding all reaction mixture to a low binding black plate**

35  $\mu$ l of HDAC assay buffer (BPS catalog number 50031)  
5  $\mu$ l of 1 mg/ml BSA  
5  $\mu$ l of 200  $\mu$ M substrate (BPS catalog number: 50037)  
5  $\mu$ l HDAC2 (5 ng/ $\mu$ l)

Always add HDAC at the last.

Incubate at 37 °C for **30 min**.

### **Setp2: stop the reaction**

add 50  $\mu$ l of fluor de lys 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)