



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

### **HDAC10**

human, recombinant, N-terminal GST Tag  
Catalog #: 50010

**Fomulated in:** 25 mM Tris-HCl, pH 8.0, 100 mM NaCl, 0.05% Tween-20 and 50% glycerol.

**Stability:** >6 months at -80°C

#### **References:**

1. Chakraborty S. (2006) J.Biol.Chem. 281(46):35070-80.
2. Wolfgang, F. (2001) J. Biol. Chem. 276(38): 35826-35835.

#### **Description:**

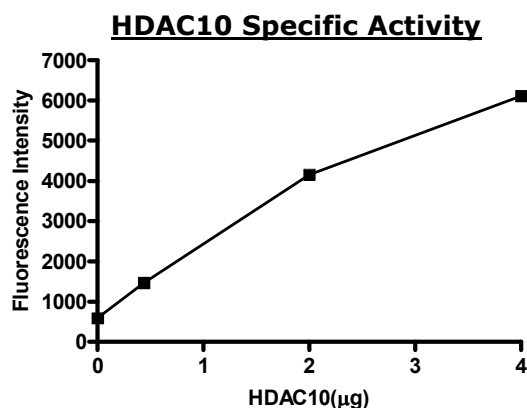
Human HDAC10 (a.a. 1-481) (GenBank Accession No. NM\_032019) with N-terminal GST tag, MW= 79 kDa, expressed in baculovirus expression system.

**Specific Activity:** 1.6 pmol/min/ug. 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 uM BPS HDAC substrate (Catalog number 50037), and 20 ng/ul HDAC10. Incubation condition: 30 min at 37°C.

#### **Application:**

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

### **Quality Assurance**

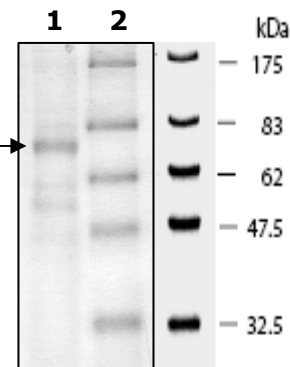


#### **10% SDS-PAGE Coomassie staining**

**Lane 1:**  
4 µg HDAC10 →

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

**MW:** 79 kDa.  
**Purity:** 70%



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 black plate**

15 µl of HDAC assay buffer (BPS catalog number 50031)  
5 µl of 1 mg/ml BSA  
5 µl of 200 uM substrate (BPS catalog number: 50037)  
25 µl of HDAC10 (0.1 µg/µl)

Always add HDAC10 enzyme at the last.

Incubate at 37 °C for 30 min.

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

add 50 µl 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)