



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 9

human, recombinant, C-terminal HIS tag  
Catalog #: 50009

Formulated in: 25 mM Tris-HCl, pH 8.0, 138 mM NaCl, 0.05% Tween-20 and 50% glycerol.

Stability: >6 months at -80°C

References:

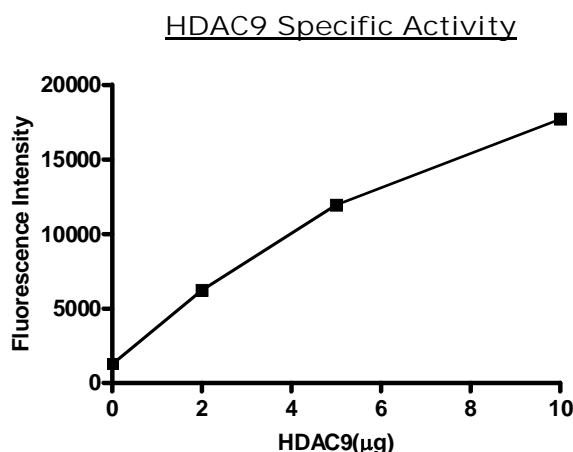
1. Han et al., (2005) J.Mol. Biol. 345, 91-102
2. Lomonte et al., (2004) J. Virol 78, 6744-6757.

Description: Human HDAC9 (GenBank Accession No. NM\_178423), amino acids 604- 1066 with C-terminal His tag, MW= 50.7 kDa, expressed in baculovirus expression system.

Specific Activity: >1000 U/ug. One U =1 pmol/min, 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 50040), and 0.2 ng/ul HDAC9. Incubation condition: 30 min at 37°C.

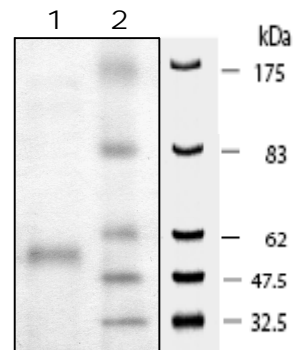
Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

### Quality Assurance



10% SDS-PAGE  
Coomassie staining

Lane 1:  
3 µg HDAC 9  
Lane 2:  
Protein Marker  
BioLabs (#P7708L)



MW: 50.7 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 50037)

### **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: 50040)  
5  $\mu$ l of HDAC9 (1 ng/ $\mu$ l)

Always add HDAC9 enzyme 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)