

Data Sheet

HDAC 9

Human, recombinant, C-terminal His-tag

Catalog #: 50009

Lot#: 130502-66 **Conc.:** 0.68 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 66 mM Imidazole, and 20% glycerol.

Stability: >6 months at -70°C . Avoid freeze/thaw cycles.

References:

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

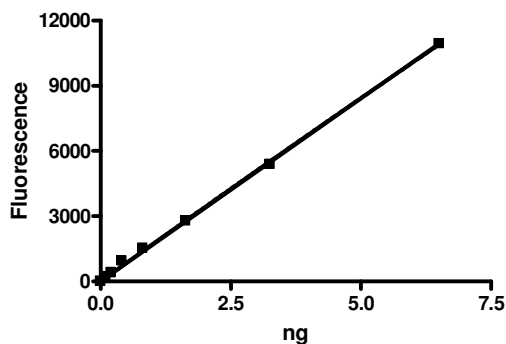
Description: Human HDAC9 (GenBank Accession No. NM_178423), amino acids 604- 1066 with C-terminal His-tag, MW= 50.8 kDa, expressed in baculovirus expression system.

Specific Activity: 3,000 pmol/min/ μg . Assay condition: 25 mM Tris/HCl, pH 8.0, 137 mM NaCl, 2.7 mM KCl, 1 mM MgCl_2 , and 0.1 mg/ml BSA, 20 μM BPS HDAC substrate (Catalog #50040), and HDAC9. 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



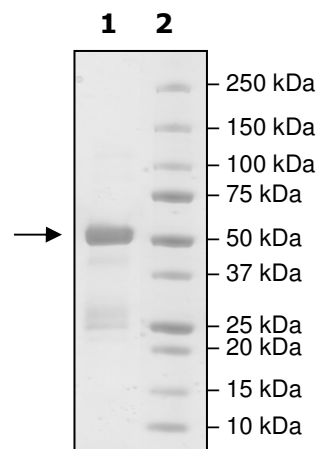
4-20% SDS-PAGE
Coomassie staining

Lane 1:
3.15 μg HDAC9

Lane 2:
Protein Marker

MW: 50.8 kDa

Purity: $\geq 77\%$



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

Please visit our website at: www.bpsbioscience.com



6044 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Sequence:

MRLVSRTHSSPAASVLPHPAMDRPLQPGSATGIAYDPLMLKHQCVCNSTTHPEHA
GRIQSIWSRLQETGLLNKCEIQGRKASLEEIQLVHSEHHSLLYGTNPLDGQKLDPRIL
LGDDSQKFFSSLPCGGLGVDSDTIWNELHSSGAARMAVGCVIELASKVASGELKNGF
AVVRPPGHHAEESTAMGFCFFNSVAITAKYLRDQLNISKILIVDLVDVHHGNGTQQAFY
ADPSILYISLHRYDEGNFFPGSGAPNEVGTGLGEGYNINIAWTGGLDPPMGDVEYLEA
FRTIVKPVAKFDPDMVLVSAGFDALEGHTPPLGGYKVTAKCFGHLTKQLMTLADGR
VVLALEGGHDLTAICDASEACVNALLGNELEPLAEDILHQSPNMNAVISLQKIIQSKY
WKSVRMVAVPRGCALAGAQLQEETETVSALASLTVDVEQPFAQEDSRTAGEPMEEE
PALHHHHHH

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
Please visit our website at: www.bpsbioscience.com