

Data Sheet

HDAC 8

Human, recombinant, C-terminal His tag

Catalog #: 50008

Lot#: 131120-1 **Conc.:** 1.1 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 200mM Imidazole, and 20% glycerol.

Stability: >6 months at -80°C

References:

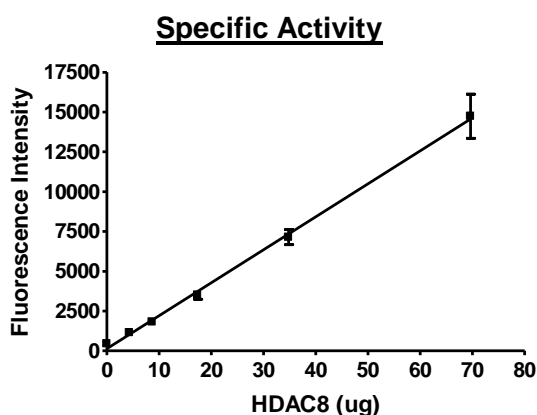
1. Lee, H. *et al.*, *Mol. Cell. Biol.* **26 (14)**, 5259-5269 (2006)
2. Gantt, S.L. *et al.*, *Biochemistry* **45 (19)**, 6170-6178 (2006)

Description: Human HDAC8 (GenBank Accession No. NM_018486), full length with C-terminal His tag, MW= 46.4 kDa, expressed in a baculovirus expression system.

Specific Activity: 300 pmol/min/μg.
 Assay condition: 25mM Tris-HCl, pH 8.0, 137mM NaCl, 2.7mM KCl, 1mM MgCl₂, 0.05% Tween 20 and 20μM HDAC class 2a substrate (50040). Incubate for 30 minutes at 37°C, followed by incubating with developer at room temperature for 20 minutes. Fluorescence intensity is measured at exc360/em460.

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

Quality Assurance

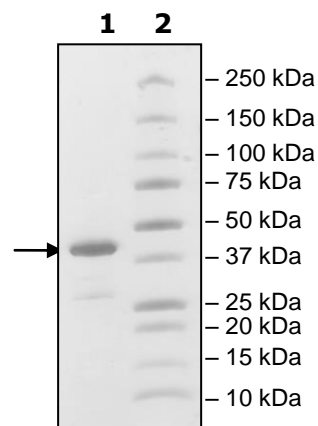


4-20% SDS-PAGE
Coomassie staining

Lane 1:
2.9 μg HDAC8

Lane 2:
Protein Marker

MW: 46.4 kDa
Purity: ≥90%



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

Amino acid sequence

NP_060956

MEEPEEPADSGQSLVPVYIYSPEYVSMCDLAKIPKRASMVHSLIEAYALHKQM
RIVKPKVASMEEMATFHTDAYLQHLQKVSQEGDDHPDSIEYGLGYDCPATEG
IFDYAAAIGGATITAAQCLIDGMCKVAINWSSGGWHHAKKDEASGFCYLNDVAVLG
ILRLRRKFERILYVDLDLHHGDGVEDAFSFTSKVMTVSLHKFSPGFFPGTGDVS
DVGLGKGRYYSVNVPIQDGIQDEKYYQICESVLKEVYQAFNPKAVVLQLGADTI
AGDPMCSFNMTVPVGIGKCLKYILQWQLATLILGGGGYNLANTARCWTYLTGVIL
GKTLSSSEIPDHEFFTAYGPDYVLEITPSCRPRDRNEPHRIQQILNYIKGNLKHVVH
HHHHH

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