



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

Data Sheet

HDAC 8

human, recombinant, C-terminal His tag
Catalog #: 50008

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:

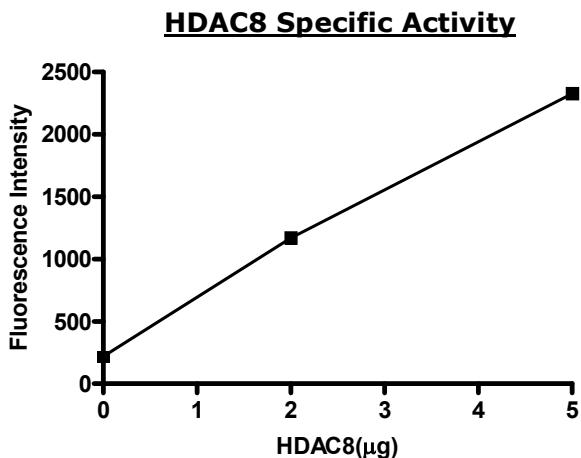
1. Lee, H. et al Mol. Cell. Biol. 26 (14), 5259-5269 (2006)
2. Gantt, S.L. et al., Biochemistry 45 (19), 6170-6178 (2006)

Description: Human HDAC8 (GenBank Accession No. NM_018486), full length with C-terminal His tag, MW= 42.4 kDa, expressed in a baculovirus expression system.

Specific Activity: 10 U/μg. One U = 1 pmol/min, Assay condition: 25 mM Tris/Cl, pH 8.0, 137 mM NaCl, 2.7 mM KCl, 1 mM MgCl₂, and 0.1 mg/ml BSA, 20 μM BPS HDAC substrate (Catalog number 50037), and 20 ng/μl HDAC8. Incubation condition: 30 min at 37°C.

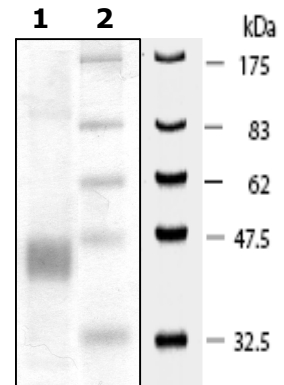
Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance



10% SDS-PAGE Coomassie staining

Lane 1:
6 μg HDAC 8
Lane 2:
Protein Marker
BioLabs (#P7708L)



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 **1.858.202.1401** or **1.858.829.3082**, Fax **1.858.259.5734**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: 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

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 μ l of HDAC assay buffer (BPS catalog number 50031)
5 μ l of 1 mg/ml BSA
5 μ l of 200 μ M substrate (BPS catalog number: 50037)
5 μ l of HDAC8 (0.2 μ g/ μ l)

Always add HDAC at the last.

Incubate at 37 °C for 30 min.

Setp2: stop the reaction

add 50 μ 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

Please visit our website at: 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

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
Please visit our website at: www.bpsbioscience.com