



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

HDAC4

human, recombinant, N-terminal GST Tag
Catalog #: 50004

Formulated 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:

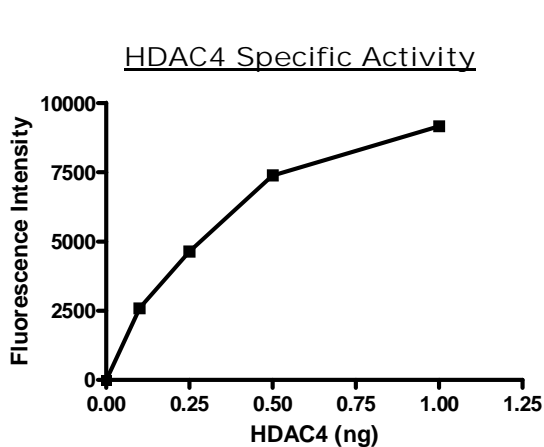
1. Leong, H. et al (2005) Mol. Endocrinol. 19, 2930-2942.
2. Zhao et al., (2005) Mol. Cell. Biol. 25, 8456-8464.

Description: Human HDAC4 (GenBank Accession No. NM_006037), amino acids 627- 1085 with N-terminal GST tag, MW= 75.2 kDa, expressed in baculovirus expression system.

Specific Activity: 14,610 pmol/min/μg, Assay condition: 25 mM Tris/Cl, pH8.0, 137 mM NaCl, 2.7 mM KCl, 1 mM MgCl₂, and 0.1 mg/ml BSA, 20 μM BPS HDAC substrate (Catalog number 50040), and 0.1 ng/μl HDAC4. Incubation condition: 30 min at 37°C.

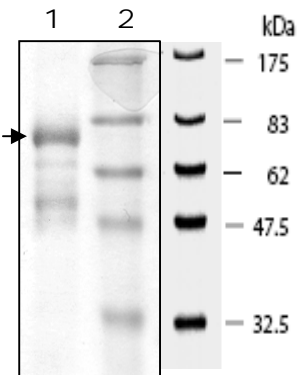
Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance



10% SDS-PAGE
Coomassie staining

Lane 1:
5 μg HDAC4 →
Lane 2:
Protein Marker
BioLabs (#P7708L)
MW: 75.2 kDa.
Purity: >50%



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

HDAC4 Fluorimetric Assay

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 ul of HDAC assay buffer (BPS catalog number 50031)
5 ul of 1 mg/ml BSA
5 ul of 200 uM substrate (BPS catalog number: 50040)
5 ul of HDAC4 (0.1 ng/ul)

Always add HDAC4 at the last.

Incubate at 37 °C for 30 min.

Setp2: stop the reaction

add 50 ul 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

Please visit our website at: www.bpsbioscience.com