



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 5 (human, recombinant, C-terminal HIS Tag)
Catalog #: 50005

Description: Human HDAC5 (GenBank Accession No. NM_001015053), amino acids 657-1123 with C-terminal His tag, MW= 51 kDa, expressed in baculovirus expression system.

Source: Sf9 cells

Formulated in: 50 mM Tris-HCl, pH 8.0, 138 mM NaCl, 10% glycerol.

Stability: >6 months at -80°C

Purity: 50% by SDS-PAGE

Specific Activity: 0.5 U/μg. One U =1 pmol/min, 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 Biomol substrate (Catalog number 50037), and 40 ng/μl HDAC5. Incubation condition: 30 min at 37°C.

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

References :

1. Spiegelberg, B.D. and Hamm, H.E. (2005) J. Biol. Chem. 280, 41769-41776
2. Waltregny et al., (2004) J. Neural Transm. 48, 273-290.

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