

Data Sheet

HDAC 5 Catalytic Domain

human, recombinant, C-terminal His-tag

Catalog #: 50005

Lot#: 100414 **Conc.:** 1.65 mg/ml

Formulated in: 45 mM Tris-HCl, pH 8.0, 124 mM NaCl, 2.4 mM KCl, 225 mM imidazole, and 10% glycerol.

Stability: >6 months at -80°C. Avoid freeze/thaw cycles.

References:

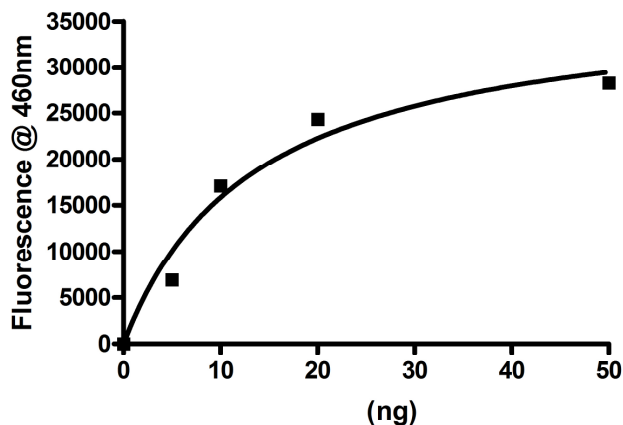
1. Spiegelberg, B.D. and Hamm, H.E. (2005) *J. Biol. Chem.* **280**, 41769-41776
2. Waltregny *et al.*, (2004) *J. Neural Transm.* **48**. 273-290.

Description: Human HDAC5 catalytic domain, variant 1 (GenBank Accession No. NM_005474), amino acid 656-1122 (end) with C-terminal His-tag, MW= 51 kDa, expressed in baculovirus expression system.

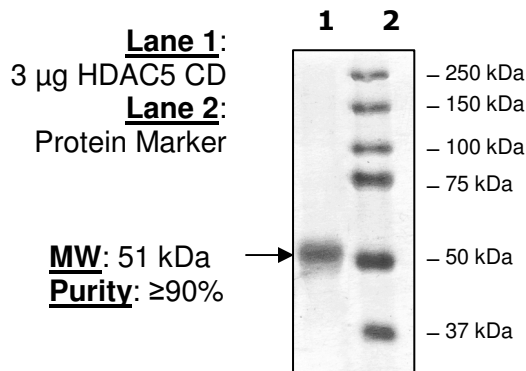
Specific Activity: 2521 pmol/min/μg. Assay condition: 25 mM Tris/HCl, pH 8.0, 137 mM NaCl, 2.7 mM KCl, 1 mM MgCl₂, and 0.1 mg/ml BSA, 30 μM BPS substrate (Catalog #50040), and HDAC5. Incubation condition: 30 min at 30°C.

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

Quality Assurance



10% SDS-PAGE Coomassie staining



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