

Data Sheet

HDAC4

Human, recombinant, N-terminal GST Tag
Catalog #: 50004
Lot#: 130115 **Conc.:** 1.5 mg/ml

Formulated in: 45 mM Tris-HCl, pH 8.0,
124 mM NaCl, 2.4 mM KCl, 18 mM
glutathione, and 10% 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.

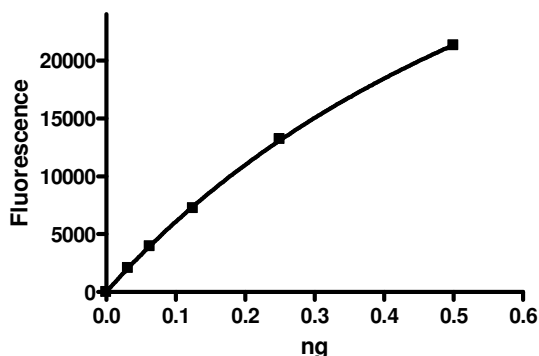
Description: Human HDAC4 (GenBank Accession No. NM_006037), amino acids 627- 1085 with N-terminal GST tag, MW= 75.2 kDa, expressed in baculovirus expression system.

Specific Activity: 103,255 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, 20 μM BPS HDAC class 2a substrate (Catalog #50040), and HDAC4. Incubation condition: 30 min at 37°C.

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

Quality Assurance

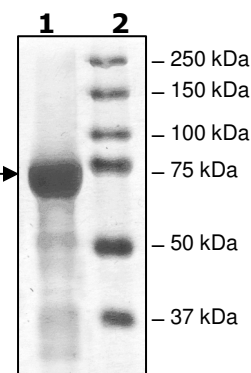
Specific Activity



10% SDS-PAGE Coomassie staining

Lane 1:
5 μg HDAC4
Lane 2:
Protein Marker

MW: 75.2 kDa
Purity: ≥60%



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