

Data Sheet

HDAC2

Human, recombinant, C-terminal FLAG-tag

Catalog #: 50052

Lot#: 140327-1

Conc.: 1.1mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, and 80 µg/ml FLAG peptide

Stability: >6 months at -80°C. Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

References:

1. Yang, W.M. *et al.*, *Proc. Natl. Acad. Sci. U.S.A.* **93 (23)**, 12845-12850 (1996).
2. Mungall, A.J. *et al.*, *Nature* **425 (6960)**, 805-811 (2003).

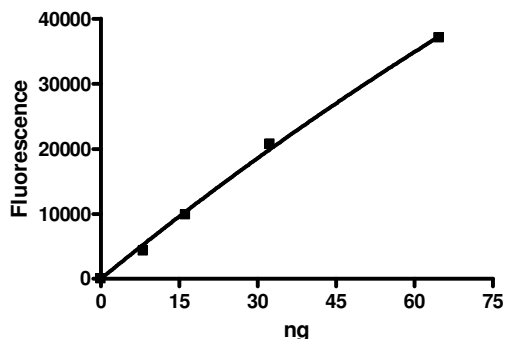
Description: Human HDAC2 (GenBank Accession No. Q92769), full length with C-terminal FLAG-tag, MW= 60 kDa, expressed in baculovirus expression system.

Specific Activity: 1020 pmol/min/µg. 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 substrate 3 (Catalog #50037), and HDAC2 (3.1 – 200 ng). Incubation condition: 30 min at 37°C followed by developing for 15 min at room temperature.

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

Quality Assurance

Specific Activity

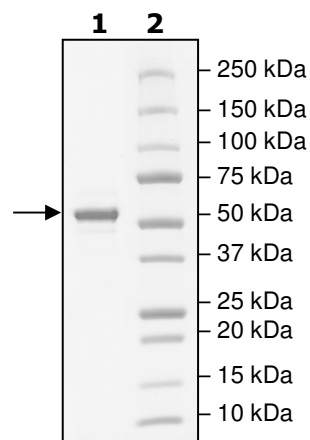


**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
2.9 µg HDAC2

Lane 2:
Protein Marker

MW: 60 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 **1.858.829.3082** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com