

Data Sheet

HDAC 7

Human, recombinant, N-terminal GST tag

Catalog #: 50007

Lot #: 150714-G2 **Conc.:** 0.06 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 250 mM NaCl, and 20% glycerol.

Stability: At 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. Chakraborty, S. (2006) *J. Biol. Chem.* **281(46)**:35070-80.
2. Wolfgang, F. (2001) *J. Biol. Chem.*

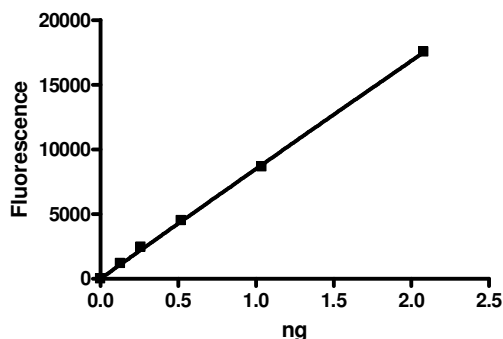
Description: Human HDAC7 (GenBank Accession No. AY302468), (a.a. 518-end) with N-terminal GST tag, MW= 78 kDa, expressed in baculovirus expression system.

Specific Activity: 16,600 pmol/min/ μg . Assay condition: 25 mM Tris/HCl, pH 8.0, 137 mM NaCl, 2.7 mM KCl, 1 mM MgCl_2 , and 0.1 mg/ml BSA, 20 μM BPS HDAC substrate 1 (Catalog #50040), and various amounts of HDAC7. Incubation conditions: 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

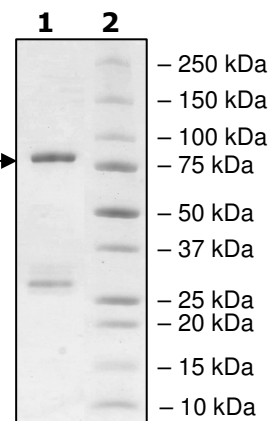
Specific Activity



4-20% SDS-PAGE Coomassie staining

Lane 1: 2 μg HDAC7
Lane 2: Protein Marker

MW: 78 kDa
Purity: $\geq 70\%$



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