

Data Sheet

HDAC10, FLAG-Tag Recombinant

Human, Recombinant, N-terminal FLAG-Tag

Catalog #: 50060

Lot #: 191121

Conc.: 0.30 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0. 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% Glycerol and 90 µg/ml FLAG peptide.

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

References:

1. Chakraborty S. (2006) *J. Biol. Chem.* **281(46)**:35070-80.

2. Wolfgang, F. (2001) *J. Biol. Chem.* **276(38)**: 35826-35835.

Description: Human HDAC10, also known as Histone Deacetylase 10, Polyamine Deacetylase HDAC10, and HD10, GenBank Accession No. NM_032019, a.a. 2-631 with N-terminal FLAG tag, expressed in a Baculovirus-infected Sf9 cell expression system. MW= 67 kDa.

Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:

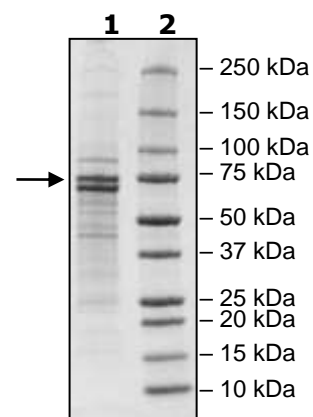
3µg HDAC10

Lane 2:

Protein Marker

MW: 67 kDa

Purity: 62%



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** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com