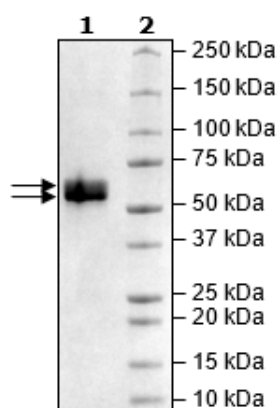


Product Information

Description:	Human recombinant HDAC1 (Homo sapiens histone deacetylase 1), full length, encompassing amino acids 1-482(end). This construct contains a C-terminal His-tag (6xHis) followed by a Flag-tag. This recombinant protein was affinity purified as a complex with endogenous tubulin.
Species	Human
Construct:	HDAC1/Tubulin (1-482(end)-His-FLAG)
Concentration:	0.84 mg/ml
Expression System:	Sf9
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 100 ng/μl FLAG peptide, and 20% Glycerol
MW:	56 kDa + Tubulin (55kDa)
Genbank Accession:	NM_004964
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
Specific Activity:	460 pmol/min/μg
Assay Conditions:	Assay was done according to HDAC1 Fluorogenic Assay Kit (BPS Bioscience #50061) with HDAC1 titrated at various concentrations.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



HDAC1 Activity

