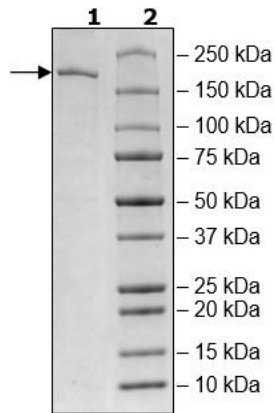


Product Information

Description:	Recombinant human HDAC6 (histone deacetylase 6), full length encompassing amino acids 1-1215(end). This protein contains an N-terminal GST-tag. A short peptide linker placed between GST and the protein includes a Thrombin Cleavage Site. This recombinant protein was affinity purified.
Background	HDAC6 is a distinct member of the histone deacetylase family, primarily found in the cytoplasm. Unlike other HDACs it targets non-histone proteins. HDAC6 regulates cell motility by deacetylating α -tubulin and helps degrade misfolded proteins through autophagy. Additionally, it influences stress responses and immune signaling by modulating acetylation of key regulatory proteins. Due to its versatile roles, HDAC6 is a promising drug target with potential therapeutic applications in cancer, neurodegenerative diseases, and immune disorders.
Species:	Human
Construct:	HDAC6 (GST-Th-1-1215(end))
Concentration:	1.05 mg/ml
Expression System:	Sf9
Purity:	$\geq 90\%$
Format:	Aqueous buffer solution.
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, and 3 mM DTT
MW:	161 kDa
Genbank Accession:	NM_006044
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:	290 pmol/min/ μg
Assay Conditions:	Assay run according to HDAC6 Fluorogenic Assay Kit (BPS Bioscience #50076) with HDAC6 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



HDAC6 Activity

