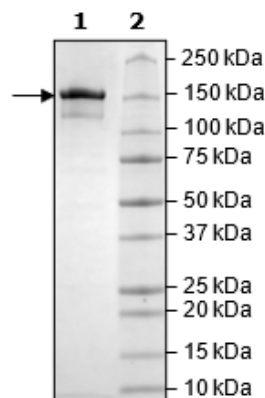


Product Information

Description:	Recombinant rat full length HDAC6 (Histone Deacetylase 6) encompassing amino acids 1-1152(end). The construct contains a N- terminal GST-Tag. This protein was affinity purified.
Background:	Histone Deacetylases (HDACs) are involved in histone deacetylation, which regulates transcription. HDAC6 is involved in multiple cellular processes, such as leptin sensitivity, cell motility, mitosis and stress granule formation. Overexpression of HDAC6 is linked to tumor formation, and mutations have been associated with Alzheimer's disease. The use of HDAC6 inhibitors presents high therapeutic potential for the treatment of neurodegenerative diseases and cancer.
Species:	Rat
Construct:	HDAC6 (GST-1-1152(end))
Concentration:	0.53 mg/ml
Expression System:	Sf9
Purity:	82%
Format:	Aqueous buffer solution.
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20 and 20% glycerol
MW:	156 kDa
Genbank Accession:	XM_003752010.4
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:	530 pmol/min/μg
Assay Conditions:	Assay was done according to HDAC6 Fluorogenic Assay Kit (BPS Bioscience #50076) with various amounts of HDAC6.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



HDAC6 Activity

