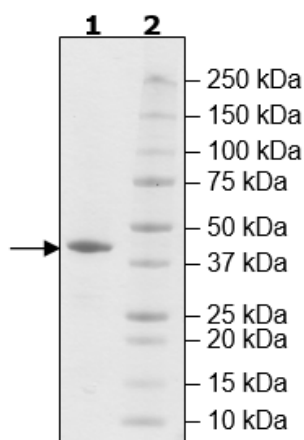


Product Information

Description:	Human histone deacetylase 8 (HDAC8), full length encompassing amino acids 1-377(end). This construct contains a C-terminal His-tag. The recombinant protein was affinity purified.
Species	Human
Construct:	HDAC8 (1-377(end)-His)
Concentration:	3.70 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, 200 mM Imidazole, and 20% glycerol
MW:	43 kDa
Genbank Accession:	NM_018486
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:	300 pmol/min/μg
Assay Conditions:	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 Fluorogenic HDAC class 2a substrate (BPS Bioscience #50040), and HDAC8 (3.1 – 200 ng). Incubation condition: 30 min at 37°C followed by developing for 15 min at room temperature.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



HDAC8 Activity

