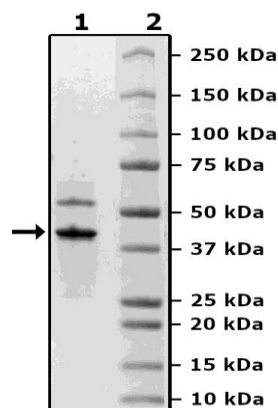


Product Information

Construct:	HDAC11 (1-347(end))
Concentration:	0.20 mg/ml
Species:	Human
Formulated In:	50 mM Tris-HCl, pH 8.0, 500 mM NaCl, 50 mM KCl, 2 mM EDTA, 3 mM DTT, 0.5% Triton-X-100, and 10% Glycerol
Expression System:	Sf9
Format:	Aqueous buffer solution.
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_024827
MW:	40 kDa
Purity:	80%
Specific Activity:	25 pmol/min/g
Assay Conditions:	Assay was done according to HDAC11 Fluorogenic Assay Kit (BPS Bioscience #50687) with various amounts of HDAC11.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



HDAC11 Activity

