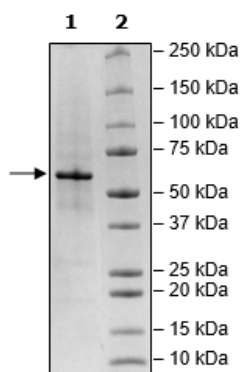


## Product Information

<b>Construct:</b>	HDAC2 (1-488(end)-His)
<b>Concentration:</b>	0.63 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol
<b>Expression System:</b>	Sf9
<b>Format:</b>	Aqueous buffer solution
<b>Stability:</b>	At least 6 months at -80°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>Genbank Accession:</b>	NM_001527
<b>MW:</b>	56 kDa
<b>Purity:</b>	85%
<b>Specific Activity:</b>	675 pmol/min/μg
<b>Assay Conditions:</b>	Assay was done according to the HDAC2 Fluorogenic Assay Kit (BPS Bioscience #50062) with HDAC2 titrated from 20 ng-0.30- ng/rxn.
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

### 4-20% SDS-Page Coomassie Staining



### Specific Activity

