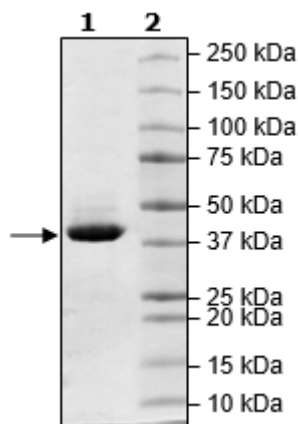


Product Information

Construct:	IDH1 (1-414(end)-FLAG)
Concentration:	2.0 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 100 µg/ml FLAG peptide, 3 mM DTT and 20% 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_005896
MW:	48 kDa
Purity:	≥90%
Specific Activity:	3,900 pmol/min/µg
Assay Conditions:	50 mM Tris (pH7.4), 10 mM MgCl ₂ , 200 µM DL-isocitrate, 200 µM NADP. Read Abs340 using Tecan Infinite 1000 (5 flash/10 ms settle time) in kinetic mode for 20 min. Molar extinction coefficient of NADPH is 6,220 M ⁻¹ cm ⁻¹
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Specific Activity

