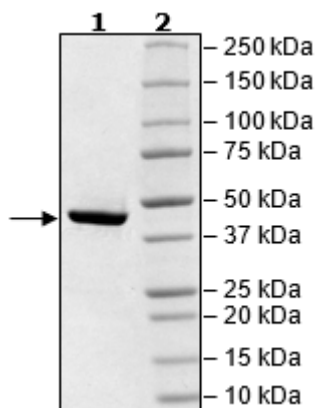


## Product Information

<b>Construct:</b>	IDH2 (R140Q) (1-452(end)-FLAG)
<b>Mutation:</b>	R140Q
<b>Concentration:</b>	0.36 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, 100 µg/ml FLAG peptide
<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_002168
<b>MW:</b>	52 kDa
<b>Purity:</b>	≥90%
<b>Specific Activity:</b>	61 pmol/min/µg
<b>Assay Conditions:</b>	Assay run according to IDH2(R140Q) Assay Kit (BPS Bioscience #79309) with IDH2 titrated at various concentrations.
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

4-20% SDS-Page Coomassie Staining



IDH2 Activity

