

IDH2 (His)

CATALOG NO.: IDH-11-315

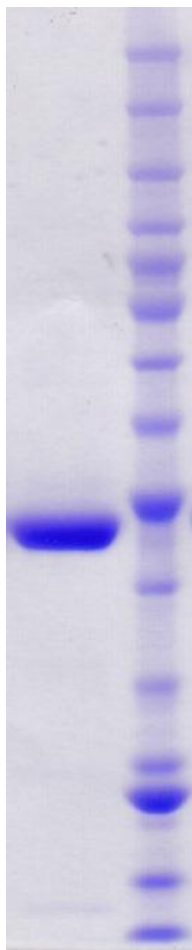
LOT NO.: 2043

DESCRIPTION: Human recombinant IDH2 (residues 40-452; Genbank Accession # NM_002168.3; MW = 47.8 kDa) expressed with a C-terminal His-tag in *E. coli*.

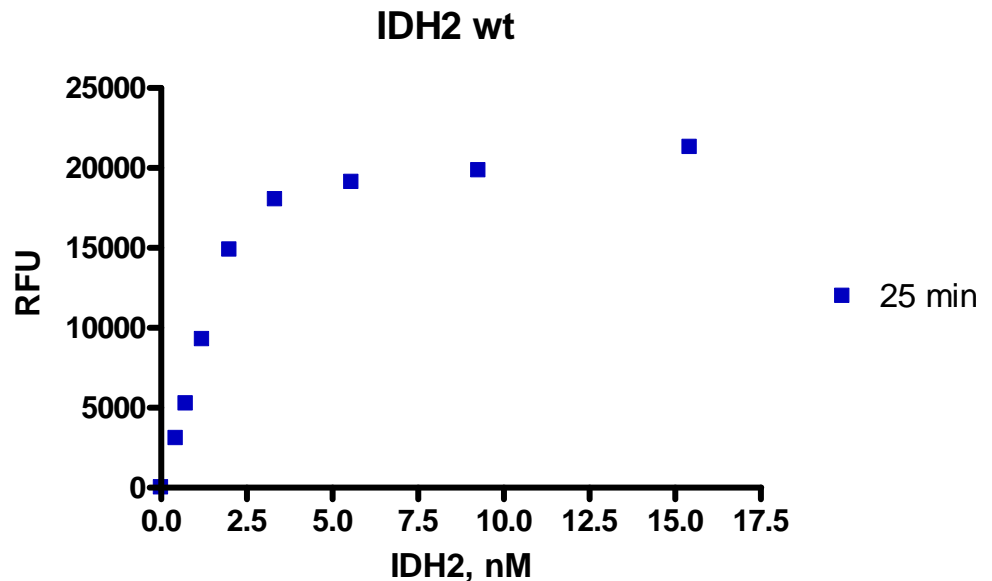
PURITY: >90% by SDS-PAGE

SUPPLIED AS: 1.44 µg/µL in 50 mM Tris HCl, pH 7.5, 500 mM NaCl, 1 mM TCEP, 10% glycerol

STORAGE: -70°C. Thaw quickly and store on ice before use. The remaining, unused, undiluted protein should be snap frozen, for example in a dry ice ethanol bath or liquid nitrogen. Minimize freeze/thaws if possible, but very low volume aliquots (<5 µl) or storage of diluted enzyme is not recommended.



Coomassie blue-stained SDS-PAGE (12% acrylamide) of 4µg of RBC IDH2 (His). MW markers (right) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, 50, 40, 30, 25, 20, 15, 10 kDa.



IDH2 wt dehydrogenase activity: NADP dependent oxidation of isocitrate was determined by quantification of produced NADPH using diaphorase/ resazurin detection. 30 uL reaction contained 25 uM NADP, 200 uM isocitrate, 15 ug/ml diaphorase, 30 uM resazurin and variable amount of IDH2 wt. After incubation at RT for 25 minutes, resulting fluorescence (528 exc/590 em) was measured using Synergy H4 plate reader (Biotek). RFU is proportional to uM of NADPH produced by enzyme.

This product is NOT intended for therapeutic or diagnostic use in animals or in humans.