

IDH2-R140K (His)

CATALOG NO.: IDH-11-353

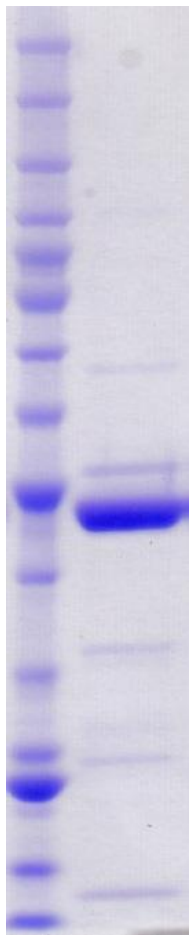
LOT NO.: 2045

DESCRIPTION: Mutant human recombinant IDH2 with lysine (K) substituted for arginine (R140) (otherwise contains wild-type residues 40-452; Genbank Accession # NM_002168.3; MW = 47.8 kDa) expressed with a C-terminal His-tag in *E. coli*.

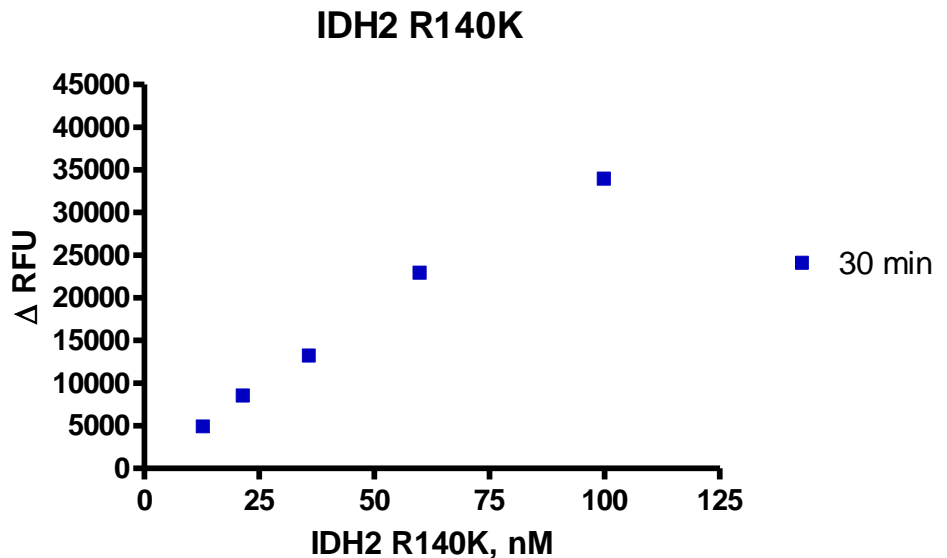
PURITY: >60% by SDS-PAGE

SUPPLIED AS: 0.77 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-R140K (His). MW markers (left) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, 50, 40, 30, 25, 20, 15, 10 kDa.



IDH2 R140K Activity Assay. NADPH-dependent reduction of α-ketoglutarate was determined by quantification of remaining NADPH using diaphorase/resazurin detection. The 20 µL reaction contained 15 µM NADPH, 10 mM α-KG and a variable amount of IDH2-R140K. After incubation at room temperature for 30 minutes, the reaction was quenched by the addition of diaphorase and resazurin (15 µg/ml and 30 µM respectively). The resulting fluorescence (ex. 528nm/em. 590nm) was measured using a Synergy H4 plate reader (Biotek). An increase in ΔRFU represents oxidation of NADPH.

Note: IDH2 R140K has wt dehydrogenase activity (not shown)

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