

IDH2-R140Q (His)

CATALOG NO.: IDH-11-351

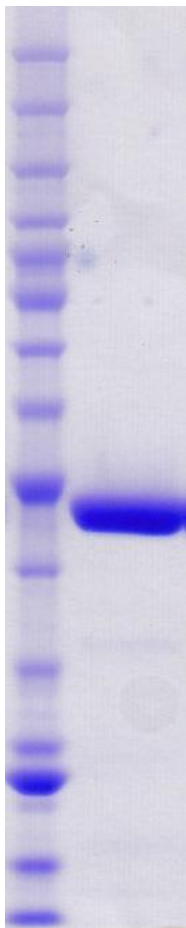
LOT NO.: 2053

DESCRIPTION: Mutant human recombinant IDH2 with glutamine (Q) 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*.

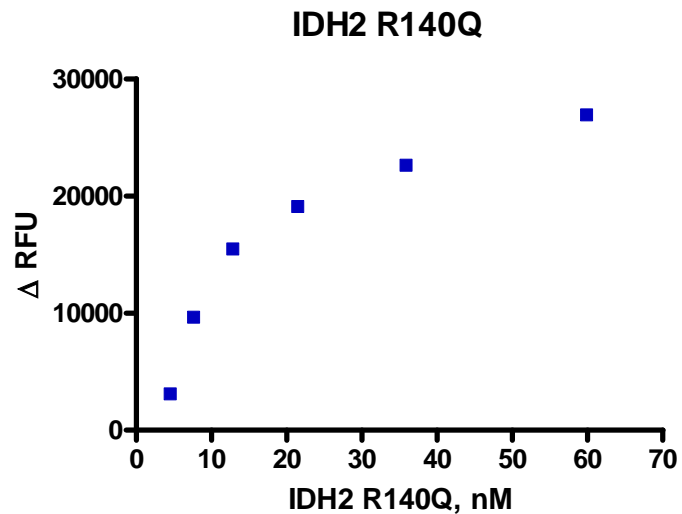
PURITY: >90% by SDS-PAGE

SUPPLIED AS: 1.86 µg/µL in 50 mM HEPES, pH 7.5, 200 mM NaCl, 3 mM DTT, 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-R140Q (His). MW markers (right) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, 50, 40, 30, 25, 20, 15, 10 kDa.



IDH2 R140Q 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-R140Q. 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, where maximum signal indicates complete oxidation of NADPH.

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