

IDH2-R172Q (His)

CATALOG NO.: IDH-11-352

LOT NO.:

DESCRIPTION: Mutant human recombinant IDH2 with glutamine (Q) substituted for arginine (R172) (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: >50% by SDS-PAGE

SUPPLIED AS: _ $\mu\text{g}/\mu\text{L}$ 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 5 μg of RBC IDH2-R172Q (His). MW markers (left) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, **50**, 40, 30, 25, **20**, 15, 10 kDa.

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