

**IDH1-R132C (His-tagged)** (Isocitrate Dehydrogenase, [NADP] cytoplasmic; Cys-132 mutant)

**CATALOG NO.:** IDH-11-324

**LOT NO.:**

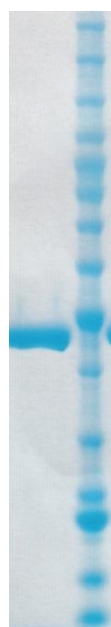
**DESCRIPTION:** Mutant human recombinant IDH1 with cysteine (C) substituted for arginine (R132) (otherwise contains wild-type residues 1-414; Genbank Accession # NM\_005896.3; MW = 47.7 kDa) expressed with a C-terminal His-tag in *E. coli*. IDH1-R132C is a cancer-associated, gain-of-function mutant form of IDH1 which can catalyze the reduction of  $\alpha$ -ketoglutarate, at the expense of NADPH oxidation, to form the 'oncometabolite' and inhibitor of various  $\alpha$ -ketoglutarate-dependent enzymes, 2-hydroxyglutarate (2-HG)<sup>1</sup>.

**PURITY:** >95% by SDS-PAGE

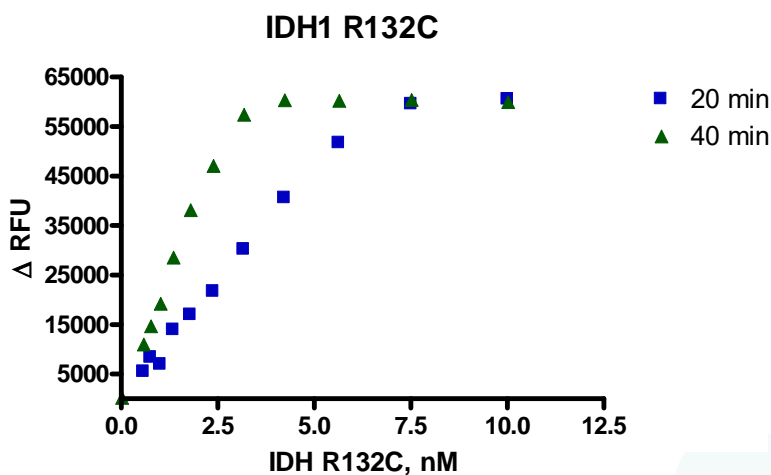
**SUPPLIED AS:**  $\_\mu\text{g}/\mu\text{L}$  in 50 mM HEPES, pH 7.5, 200 mM NaCl, 3 mM DTT, 10% glycerol

**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  $\mu\text{L}$ ) or storage of diluted enzyme is not recommended.

**REFERENCE:** 1) L. Dang *et al. Nature* 2009 **462** 739;



Coomassie blue-stained SDS-PAGE (12% acrylamide) of 2  $\mu\text{g}$  of RBC IDH1-R132C (His). MW markers (right) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, 50, 40, 30, 25, 20, 15, 10 kDa.



**IDH1 R132C Activity Assay.** NADPH-dependent reduction of  $\alpha$ -ketoglutarate was determined by quantification of remaining NADPH using diaphorase/resazurin detection. The 20  $\mu\text{L}$  reaction contained 15  $\mu\text{M}$  NADPH, 0.5 mM  $\alpha$ -KG and a variable amount of IDH1 R132C. After incubation at room temperature for 20 or 40 minutes, the reaction was quenched by the addition of diaphorase and resazurin (15  $\mu\text{g}/\text{ml}$  and 25  $\mu\text{M}$  respectively). The resulting fluorescence (ex. 528nm/em. 590nm) was measured using a Synergy H4 plate reader (Biotek). An increase in  $\Delta\text{RFU}$  represents oxidation of NADPH, where the maximum signal (~65,000  $\Delta\text{RFU}$ ) represents complete oxidation of 15  $\mu\text{M}$  NADPH.

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

Reaction Biology

1 Great Valley Parkway, Malvern PA, USA 19355

requests@reactionbiology.com www.reactionbiology.com