

PRODUCT INFORMATION



IDH1 (R132H mutant; human, recombinant)

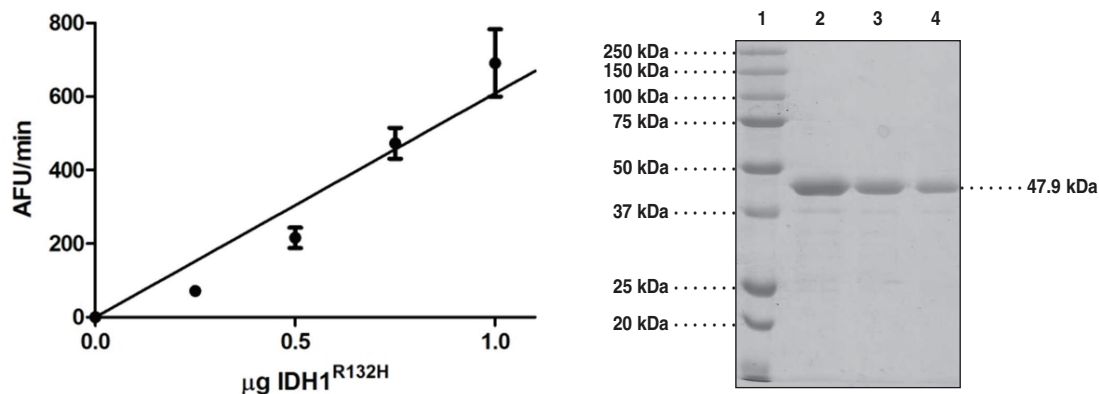
Item No. 14132

Overview and Properties

Synonym:	Isocitrate Dehydrogenase (NADP) Cytoplasmic
Source:	Active recombinant C-terminal histidine-tagged protein expressed in <i>E. coli</i>
Amino Acids:	2-414 (full-length)
Uniprot No.:	O75874
Molecular Weight:	47.9 kDa
Storage:	-80°C (as supplied)
Stability:	≥6 months
Purity:	batch specific (≥95% estimated by SDS-PAGE)
Supplied in:	50 mM Tris-HCl, pH 7.5, with 200 mM sodium chloride, 5 mM β-mercaptoethanol, and 10% glycerol
Protein	
Concentration:	batch specific mg/ml
Activity:	batch specific U/ml
Specific Activity:	batch specific U/mg
Unit Definition:	One unit is defined as the amount of enzyme required to convert 1 nmol of NADPH to NADP ⁺ , using 15 mM α-ketoglutarate as a substrate, per minute at room temperature in 25 mM Tris-HCl, pH 7.5, 150 mM sodium chloride, and 5 mM MgCl ₂ .

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Images



Lane 1: MW Markers
Lane 2: IDH1 R123H (5 μg)
Lane 3: IDH1 R123H (2.5 μg)
Lane 4: IDH1 R123H (1.25 μg)

Representative gel image shown; actual purity may vary between each batch.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 05/06/2024

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Isocitrate dehydrogenases (IDHs) are nicotinamide adenine dinucleotide (NAD⁺) and NAD phosphate (NADP⁺)-dependent enzymes that catalyze the third step of the tricarboxylic acid cycle. IDHs catalyze oxidative decarboxylation of isocitrate producing α -ketoglutarate (α -KG) and carbon dioxide. IDH1 (cytosolic) and IDH2 (mitochondrial) are NADP⁺-dependent enzymes that catalyze reversible reactions. The IDH3 isoform, a NAD⁺-dependent multisubunit enzyme, is irreversible and allosterically regulated by a variety of positive (calcium, ADP, and citrate) and negative (adenosine triphosphate, NADH, and NADPH) effectors.¹ IDH1 and IDH2 are mutated in >70% of lower grade gliomas.² The most common IDH mutation, Arg132His, imparts new gain of function catalytic activity leading to the NADPH-dependent conversion of α -KG to 2-hydroxyglutarate.^{3,4} Astrocytes expressing IDH1 R132H mutant have been shown to produce markedly increased levels of the R-2-hydroxyglutarate enantiomer, leading to transformation of cells through the hypoxia-inducible factor prolyl 4-hydroxylase, EGLN.⁵

References

1. Raimundo, N., Baysal, B.E., and Shadel, G.S. Revisiting the TCA cycle: Signaling to tumor formation. *Trends Mol. Med.* **17(11)**, 641-649 (2011).
2. Turcan, S., Rohle, D., Goenka, A., *et al.* IDH1 mutation is sufficient to establish the glioma hypermethylator phenotype. *Nature* **483(7390)**, 479-483 (2012).
3. Reitman, Z.J. and Yan, H. Isocitrate dehydrogenase 1 and 2 mutations in cancer: Alterations at a crossroads of cellular metabolism. *J. Natl. Cancer Inst.* **102(13)**, 932-941 (2010).
4. Dang, L., White, D.W., Gross, S., *et al.* Cancer-associated IDH1 mutations produce 2-hydroxyglutarate. *Nature* **462(7274)**, (2009).
5. Koivunen, P., Lee, S., Duncan, C.G., *et al.* Transformation by the (R)-enantiomer of 2-hydroxyglutarate linked to EGLN activation. *Nature* 1-7 (2012).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM