

# PRODUCT INFORMATION

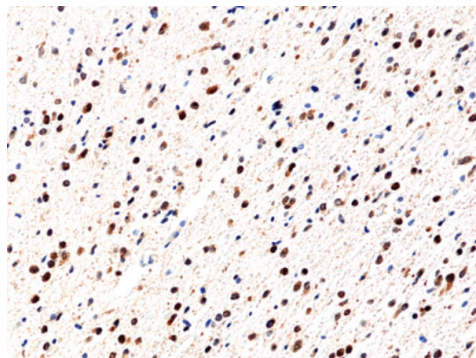
## IDH1 R132H Rabbit Monoclonal Antibody (RM390)

Item No. 32326

### Overview and Properties

<b>Contents:</b>	This vial contains 100 µl of protein A-affinity purified monoclonal antibody.
<b>Synonym:</b>	Isocitrate Dehydrogenase 1
<b>Immunogen:</b>	Peptide corresponding to the IDH1 R132H mutant
<b>Cross Reactivity:</b>	(+) IDH R132H; (-) IDH without the R132H mutation
<b>Species Reactivity:</b>	(+) Human
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥1 year
<b>Storage Buffer:</b>	PBS, with 50% glycerol, 1% BSA, and 0.09% sodium azide
<b>Clone:</b>	RM390
<b>Host:</b>	Rabbit
<b>Isotype:</b>	IgG
<b>Applications:</b>	Immunohistochemistry (IHC); the recommended starting dilution is 1:100-1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Image



Immunohistochemical staining of formalin-fixed and paraffin-embedded human glioblastoma tissue using IDH1 R132H Rabbit Monoclonal Antibody (RM390) at a dilution of 1:100.

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, 02/10/2024

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 dehydrogenase 1 (IDH1) is a NADP<sup>+</sup>-dependent enzyme and a member of the IDH family.<sup>1</sup> It is composed of a large domain, a small domain, and a clasp domain and exists as an asymmetric homodimer with active-site clefts formed by the large and small domains of one monomer and the small domain of the other. IDH1 is localized to the cytoplasm and catalyzes the oxidative decarboxylation of isocitrate to  $\alpha$ -ketoglutarate ( $\alpha$ -KG) using NADP<sup>+</sup> as a cofactor.<sup>2</sup> A mutant form of IDH1, where arginine is replaced by histidine at position 132 (IDH1<sup>R132H</sup>), changes the conformation of the IDH1<sup>R132H</sup> homodimer and IDH1<sup>R132</sup>/IDH1<sup>R132H</sup> heterodimer.<sup>3,4</sup> The IDH1<sup>R132</sup>/IDH1<sup>R132H</sup> heterodimer is unable to achieve the closed conformation needed to produce  $\alpha$ -KG and is enzymatically inactive. The IDH1<sup>R132H</sup> homodimer is not only enzymatically inactive for its native function but gains the ability to reduce  $\alpha$ -KG to D-hydroxyglutarate (D-2-HG) using NADPH as a cofactor. Exogenous IDH1<sup>R132H</sup> is oncogenic *in vitro* and the expression of IDH1<sup>R132H</sup> in the mouse brain induces tumor nodules in a process resembling gliomagenesis.<sup>5,6</sup> The IDH1<sup>R132H</sup> mutation has frequently been found in tumor tissue derived from patients with secondary glioblastomas but is associated with increased survival.<sup>5</sup> Cayman's IDH1 R132H Rabbit Monoclonal Antibody (RM390) can be used for immunohistochemistry (IHC) applications.

## References

1. Xu, X., Zhao, J., Xu, Z., *et al.* Structures of human cytosolic NADP-dependent isocitrate dehydrogenase reveal a novel self-regulatory mechanism of activity. *J. Biol. Chem.* **279**(32), 33946-33957 (2004).
2. Frezza, C., Tennant, D.A., and Gottlieb, E. IDH1 mutations in gliomas: When an enzyme loses its grip. *Cancer Cell* **17**(1), 7-9 (2010).
3. Yang, B., Zhong, C., Peng, Y., *et al.* Molecular mechanisms of "off-on switch" of activities of human IDH1 by tumor-associated mutation R132H. *Cell Res.* **20**(11), 1188-1200 (2010).
4. Zhao, S. and Guan, K.-L. IDH1 mutant structures reveal a mechanism of dominant inhibition. *Cell Res.* **20**(12), 1279-1281 (2010).
5. Huang, L.E. Friend or foe—IDH1 mutations in glioma 10 years on. *Carcinogenesis* **40**(11), 1299-1307 (2019).
6. Bardella, C., Al-Dalahmah, O., Krell, D., *et al.* Expression of Idh1<sup>R132H</sup> in the murine subventricular zone stem cell niche recapitulates features of early gliomagenesis. *Cancer Cell* **30**(4), 578-594 (2016).

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