

# IDH1 monoclonal antibody

**Monoclonal antibody to IDH1 R132H mutation for use with IHC applications**

IDH1 (Isocitrate Dehydrogenase 1) is an enzyme found in the cytoplasm and peroxisomes. IDH1 is a metabolic gene that is often mutated in human cancer. It is involved in metabolic and cellular pathways, specifically the Krebs cycle. Mutations are frequently found in diffuse low-grade gliomas and secondary glioblastomas, but rarely in primary glioblastomas. The most notable mutation in this enzyme is replacing amino acid arginine with histidine at codon 132 (R132H).

This antibody is covered by our [Worry-Free Guarantee](#).

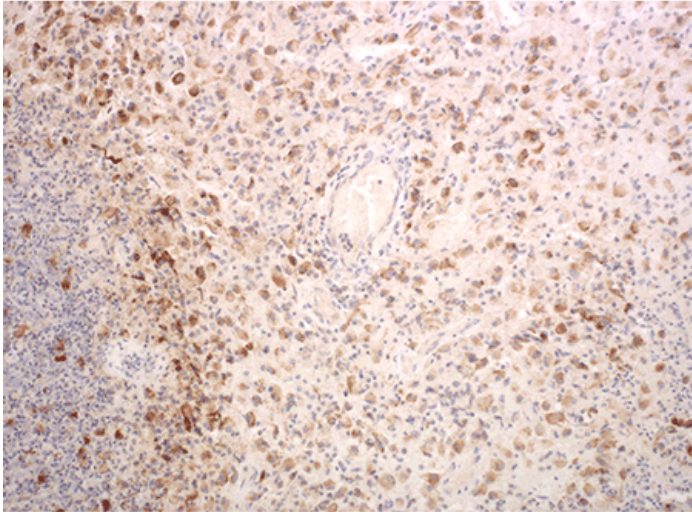
## Ordering Information

[Order Online »](#)

ENZ-ABS774-0100	100µl
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)



Formalin-fixed paraffin-embedded glioblastoma tissue stained using Enzo's IDH1 antibody (Prod. No. ENZ-ABS774). IHC staining was performed on an automated Leica BOND III.

# Handling & Storage

Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	+4°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

# Product Details

Alternative Name	Isocitrate dehydrogenase 1
Application	IHC
Application Notes	<p>The IDH1 antibody can be used to qualitatively identify the presence of IDH1 R132H protein mutation in FFPE tissue sections using IHC test methods.</p> <p>For IHC applications: Dilute the concentrate as appropriate in antibody diluent, and then apply 150 µL to the slide, as indicated in the user's IHC protocol.</p>
Formulation	Liquid. In Tris buffer, pH 7.5, containing 1% BSA and <0.1% sodium azide.
Host	Mouse
Immunogen	R132H mutant site
Isotype	IgG
Purity Detail	Affinity-purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (1:50-1:200)Optimal conditions must be determined individually for each application.
Species Reactivity	Human
UniProt ID	O75874

Worry-free Guarantee

This antibody is covered by our [Worry-Free Guarantee](#)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)