

IDH1-R132H

monoclonal antibody

(HMAb-1)

Anti-human isocitrate dehydrogenase 1 (IDH1) - R132H antibody binds to IDH1-mutated protein, but does not bind the wild-type IDH1 protein. Mutations in the IDH 1 gene are commonly found in human glioma, with the majority of low-grade gliomas harboring a recurrent point mutation IDH1-R132H. IDH1-R132H can be used as a diagnostic marker to help differentiate infiltrating gliomas from gliosis, and as a prognostic marker for gliomas and secondary glioblastoma multiforme.

Isocitrate dehydrogenase 1 (NADP+), soluble is an enzyme that in humans is encoded by the IDH1 gene. Isocitrate dehydrogenases catalyze the oxidative decarboxylation of isocitrate to 2-oxoglutarate. IDH1 is one of three isocitrate dehydrogenase isozymes, the other two being IDH2 and IDH3. IDH1 mutations are heterozygous, typically involving an amino acid substitution in the active site of the enzyme in codon 132. The mutation results in a loss of normal enzymatic function and the abnormal production of 2-hydroxyglutarate (2-HG). Mutations in IDH1 are often the first hit in the development of diffuse gliomas, suggesting IDH1 mutations as key events in the formation of these brain tumors.

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

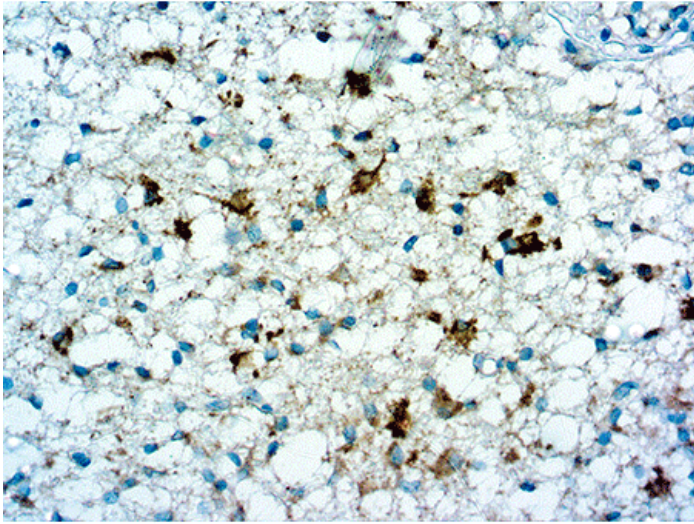
Ordering Information

[Order Online »](#)

| | |
|-----------------|-------|
| ENZ-ABS853-1000 | 1ml |
| ENZ-ABS853-0100 | 100µl |

Manuals, SDS & CofA

[View Online »](#)



Formalin fixed paraffin embedded anaplastic
oligodendroglioma stained with IDH1 antibody

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name Isocitrate dehydrogenase 1

Application IHC (PS)

Application Notes Pretreatment: Tris/EDTA Buffer pH 9.0

Clone HMAb-1

Formulation This product is supplied as a purified tissue culture supernatant and contains sodium azide as a preservative.

Host Mouse

Immunogen KLH-conjugated linear peptide corresponding to human Isocitrate Dehydrogenase 1 (IDH1).

Isotype IgG1k

Positive Control Human Glioma tissue

Recommendation Dilutions/Conditions Antibody dilutions should be appropriately adjusted and verified according to the detection system used. Optimal conditions must be determined individually for each application.

Species Reactivity Human

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

.



ENZO LIFE SCIENCES,
INC.
Phone: +1 940 948 9120
info@enzolifesciences.com
usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info@enzolifesciences.com
be@enzolifesciences.com

France
Phone: +33 472 440 655
info@enzolifesciences.com
fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info@enzolifesciences.com
de@enzolifesciences.com

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