

ALDH1L1 polyclonal antibody

The anti-ALDH1L1 antibody enables specific detection of ALDH1L1, a key astrocyte marker and enzyme involved in folate metabolism and detoxification, optimized for IHC applications in neurobiology and metabolic research.

ALDH1L1 is an enzyme in the aldehyde dehydrogenase superfamily that plays a critical role in folate metabolism by catalyzing the oxidation of 10-formyltetrahydrofolate to carbon dioxide and tetrahydrofolate. It is highly expressed in astrocytes, making it a reliable and specific marker for studying astrocytic populations in the central nervous system. ALDH1L1 is also involved in cellular detoxification processes, protecting cells from oxidative damage by metabolizing aldehydes generated during lipid peroxidation and other oxidative stress pathways.

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

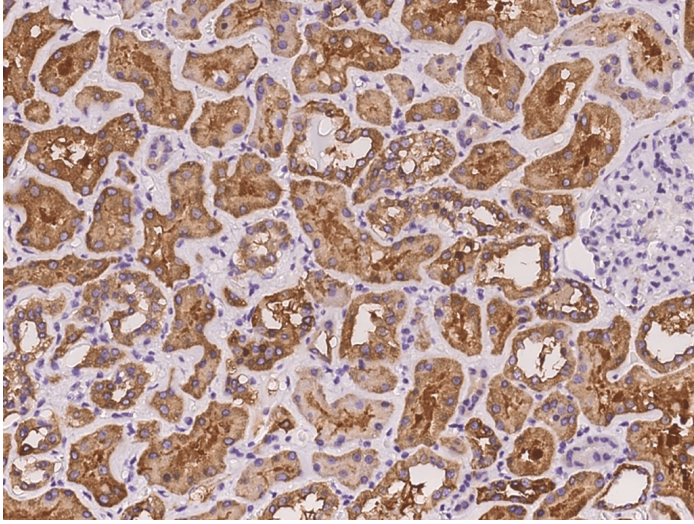
Ordering Information

[Order Online »](#)

ENZ-ABS1055-0100	100µl
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Immunohistochemical staining of human ALDH1L1 in human kidney with rabbit polyclonal antibody at 1:100 dilution, formalin-fixed paraffin embedded sections.

Handling & Storage

Handling Avoid freeze-thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Aldehyde dehydrogenase 1 family member L1, Cytosolic 10-formyltetrahydrofolate dehydrogenase

Application IHC (PS)

Formulation Liquid. 0.2 µm filtered solution in PBS. Carrier-free. Preservative-free.

Host Rabbit

Immunogen E. coli-derived human ALDH1L1 fragment.

Isotype IgG

Purity Detail Protein A and antigen affinity purified.

Recommendation Immunohistochemistry (paraffin sections) (1:50-1:200)

Dilutions/Conditions Suggested dilutions/conditions may not be available for all applications. 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: 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)