

Olig-2 recombinant monoclonal antibody (064)

The anti-Olig2 antibody is a reliable tool for detecting human Olig2, a basic helix-loop-helix transcription factor crucial for the development and maintenance of oligodendrocytes and motor neurons.

Olig-2 plays a key role in the specification and differentiation of oligodendrocytes and motor neurons in the central nervous system (CNS). Olig-2 is essential for the formation of myelin-producing oligodendrocytes, which are vital for proper neuronal function and axonal insulation. In neurodevelopment, Olig-2 regulates neural progenitor proliferation and lineage specification, while in disease contexts, it is associated with gliomas, particularly oligodendrogliomas, where it serves as a diagnostic marker. Olig-2's involvement in CNS injury and repair also highlights its importance in neurodegeneration and regenerative medicine.

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

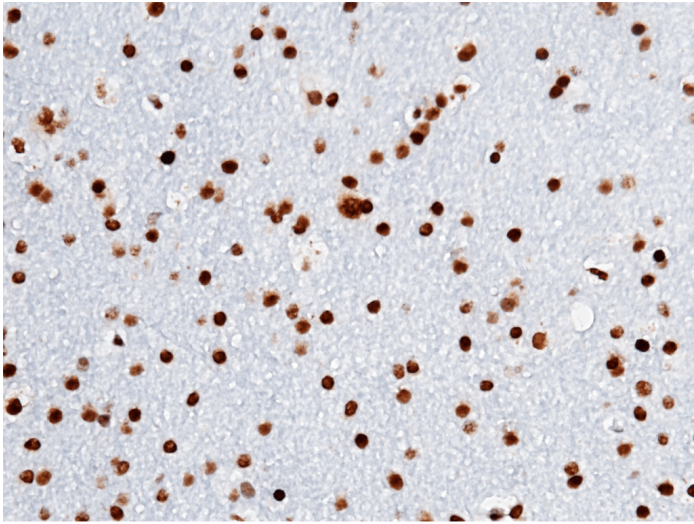
Ordering Information

[Order Online »](#)

ENZ-ABS1043-0100	100µl
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Immunohistochemical staining of human olig-2 in human brain with rabbit monoclonal antibody (1:200, 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 Oligo 2, Oligodendrocyte transcription factor 2

Application IHC (PS), WB

Clone 064

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

Host Rabbit

Immunogen Synthetic peptide corresponding to the N-terminus of the Human Olig-2.

Isotype IgG

Purity Detail Protein A purified.

Recommendation Immunohistochemistry (paraffin sections) (1:100-1:500)

Dilutions/Conditions Western Blot (1:500-1:2000)

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-
fr@enzolifesciences.com](mailto:info-fr@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)