# 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**

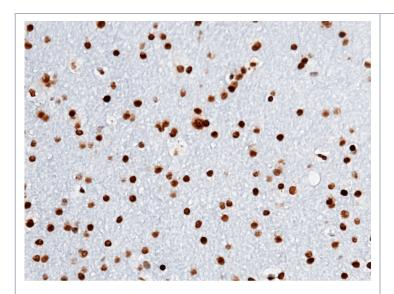
Order Online »

ENZ ADC4	042 0400	4
ENZ-ABS1	U43-U1UU	

100µl

Manuals, SDS & CofA

**View Online »** 



Immunohistochemical staining of human olig-2 in human brain with rabbit monoclonal antibody (1:200, formalinfixed 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 Dilutions/Conditions** Immunohistochemistry (paraffin sections) (1:100-1:500)

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

.

Last modified: January 9, 2025



info-

eu@enzolifesciences.com