DCX recombinant monoclonal antibody (01)

The anti-DCX antibody is designed for the detection of doublecortin (DCX), a microtubule-associated protein essential for neuronal migration and cortical development.

DCX (doublecortin) is a microtubule-associated protein critical for neuronal migration and cortical development. It is highly expressed in immature neurons during early brain development, where it facilitates cytoskeletal remodeling essential for neuronal dispersion and cortex lamination. DCX is widely used as a marker of neurogenesis and has been implicated in disorders such as lissencephaly X-linked type 1 and subcortical band heterotopia, which result from defects in neuronal migration. Its expression is largely restricted to the developing brain, making it a valuable target for studies of neurodevelopment and related pathologies.

This antibody is covered by our Worry-Free Guarantee.

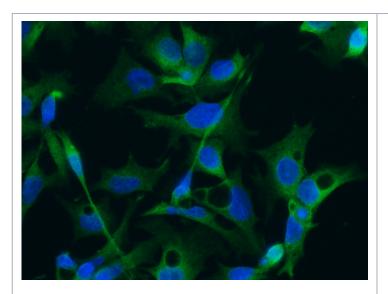
Ordering Information

Order Online »

ENZ-ABS1051-0100	100µl
ENZ-ABS1051-0100	100µ

Manuals, SDS & CofA

View Online »



Immunofluorescence staining of Human DCX in HepG2 cells. Cells were fixed with 4% PFA, permeabilized with 0.3% Triton X-100 in PBS, blocked with 10% serum, and incubated with Mouse anti-Human DCX monoclonal antibody (1:60) at 37°C 1 hour. Then cells were stained with the Alexa Fluor® 488-conjugated Goat Anti-mouse IgG secondary antibody (green) and counterstained with DAPI (blue). Positive staining was localized to cytoplasm.

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 Neuronal migration protein doublecortin

Application ICC, IF

Clone 01

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

Preservative-free.

Host Mouse

Immunogen Recombinant human doublecortin/DCX Protein.

lgG lgG

Purity Detail Protein A purified.

Recommendation Dilutions/Conditions Immunocytochemistry/Immunofluorescence (1:20-1:100)

Suggested dilutions/conditions may not be available for all

applications.

Optimal conditions must be determined individually for

each application.

Source This product is a recombinant monoclonal antibody

expressed from HEK293 cells.

Species Reactivity Human

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

.

Last modified: January 9, 2025



info-

eu@enzolifesciences.com