

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](#).

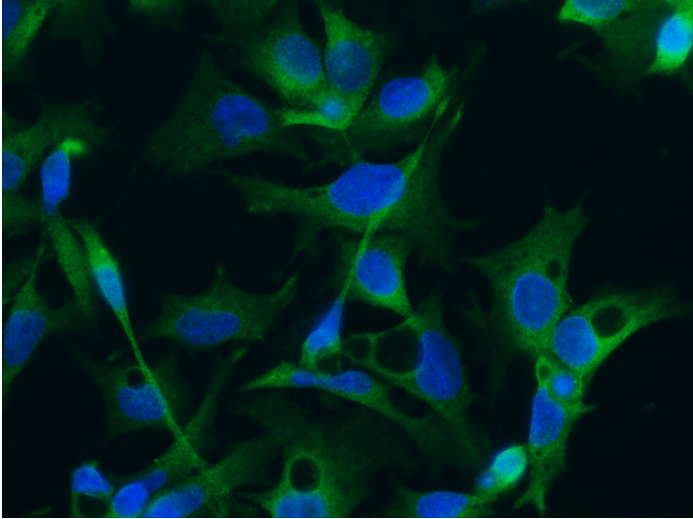
Ordering Information

[Order Online »](#)

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

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.

Isotype IgG

Purity Detail Protein A purified.

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

Dilutions/Conditions
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](#).



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)