

IDO1 (mouse) monoclonal antibody (mIDO-48)

Indoleamine 2,3-dioxygenase (IDO) is an IFN- γ inducible gene. It catalyzes the degradation of the essential amino acid L-tryptophan to N-formylkynurenine. IDO has been implicated in protection against intracellular and extracellular pathogens. The mIDO-48 monoclonal antibody recognizes mouse IDO also known as Indolamine 2,3-dioxygenase, Indole 2,3-dioxygenase, and Indoleamine-pyrrole 2,3-dioxygenase. This antibody does not appear to react with human IDO. IDO is a ubiquitously expressed cytoplasm protein with a predicted molecular weight approximately 45 kDa. Indoleamine 2,3-dioxygenase (IDO) is one the best known IFN- γ inducible genes. The product of IDO gene catalyzes the degradation of the essential amino acid L-tryptophan to N-formylkynurenine. IDO has been implicated in protection against intracellular and extracellular pathogens. It also has been shown to maintain the special immune suppressive status of immune-privileged sites such as the brain, eyes, kidney, and placenta. The IDO antibody has been shown to be useful for western blotting and Immunohistochemistry.

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

Citations: 9

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-901-0100	100 μ g
------------------	-------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	INDO, Indoleamine 2,3-dioxygenase, Indole 2,3-dioxygenase, Indoleamine-pyrrole 2,3-dioxygenase
Application	WB
Application Notes	Cited applications include IHC.
Clone	mIDO-48
Crossreactivity	Does not cross-react with IDO2 .
Formulation	Liquid. In PBS, pH 7.2, containing 0.09% sodium azide.
Gene/Protein Identifier	15930 (Entrez GeneID)
Host	Rat
Immunogen	Recombinant mouse IDO1.
Isotype	IgG2b
Purity Detail	Affinity purified.
Species Reactivity	Mouse
UniProt ID	P28776
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee



Last modified: May 29, 2024

ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430

[info-
be@enzolifesciences.com](mailto:info-be@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-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)