

AIM2 (human) monoclonal antibody (3B10)

AIM2 is a strong candidate as a tumor suppressor as it is a member of the HIN-200 (hemopoietic interferon-inducible, nuclear proteins containing a 200 amino acid repeat) family of proteins that have been shown to regulate cell growth and survival. The HIN-200 family of proteins consists of four members in human, IFI 16, MNDNA, AIM2 and IFIXI. It has been reported that AIM2 is a receptor for cytosolic dsDNA, which forms an inflammasome complex with ASC to activate caspase-1-mediated processing of IL-1 β .

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

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-891-C100	100 μ g
------------------	-------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Short Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Interferon-inducible protein AIM2, Absent in melanoma 2

Application IF, IP, WB

Application Notes Detects a band of ~39kDa by Western blot.

Clone 3B10

Formulation Liquid. In PBS, pH 7.2, containing 0.09% sodium azide.

Gene/Protein Identifier 9447 (Entrez GeneID)

Host Mouse

Immunogen Recombinant human AIM2 (aa 1-186) with GST-tag.

Isotype IgG1

Purity Detail Affinity purified.

Recommendation Western blot (0.5-2 µg/ml) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Species Reactivity Human

UniProt ID O14862

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)