

A20 monoclonal antibody (59A426)

The zinc finger protein A20 is a tumor necrosis factor and Interleukin-1 (IL-1)-inducible protein that negatively regulates NF- κ B gene expression. A20 is expressed at low levels in most organs, the highest expression levels being in lymphoid tissues. It has been shown that the A20 protein level is increased in lung neutrophils after endotoxemia.

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

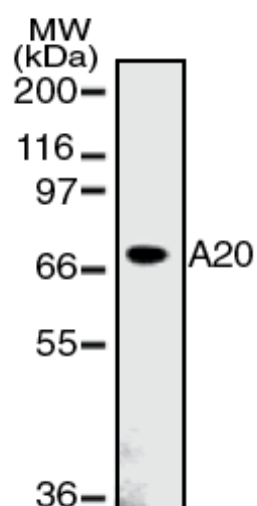
Ordering Information

[Order Online »](#)

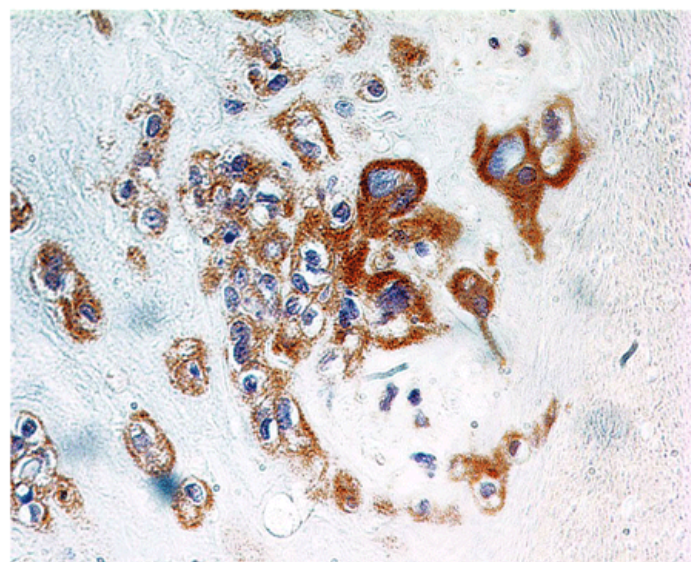
ADI-AAM-009-E	100 μ g
---------------	-------------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of human Daudi cell lysate, probed with AAM-009.



Immunohistochemistry analysis: Formalin-fixed, paraffin-embedded human placenta probed with A20, mAb (59A426) at 5 μ g/ml, cytoplasmic staining of decidual cells is observed.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application IHC (PS), IP, WB

Clone 59A426

Formulation Liquid. In PBS containing 0.05% sodium azide.

Host Mouse

Immunogen Recombinant human A20.

Isotype IgG1

Purity Detail Protein G affinity purified.

Recommendation Dilutions/Conditions Immunohistochemistry (5µg/ml)Immunoprecipitation (1-2µg/ml)Western Blot (3-5µg/ml, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Human

UniProt ID P21580

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-
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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-fr@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)