

ASC polyclonal antibody

ASC (Apoptosis-associated speck-like protein containing a CARD domain) is a bipartite protein comprising two protein-protein interaction domains, a Pyrin domain (PYD) and a caspase recruitment domain (CARD). Proteins containing these domains play pivotal roles in regulating apoptosis and immune response pathways. Reported functions of ASC include apoptosis, activation of inflammatory caspases and regulation of NFκB.

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

Citations: 39

[View Online »](#)

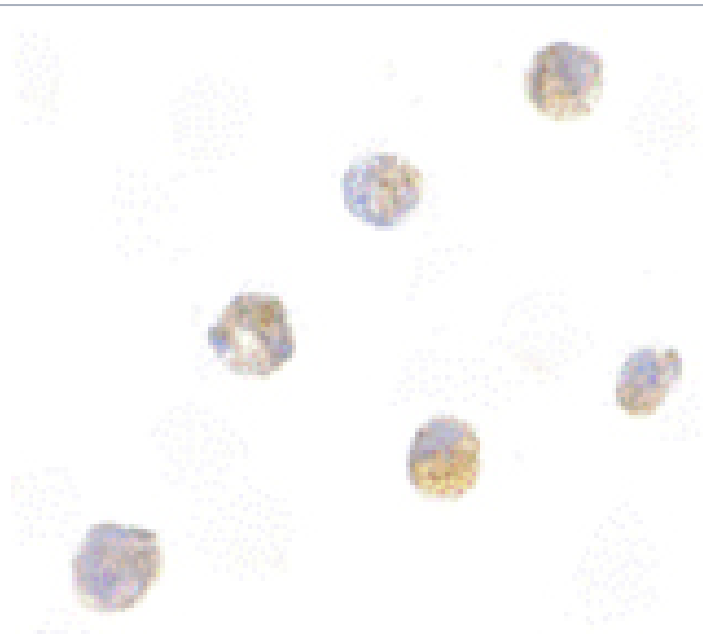
Ordering Information

[Order Online »](#)

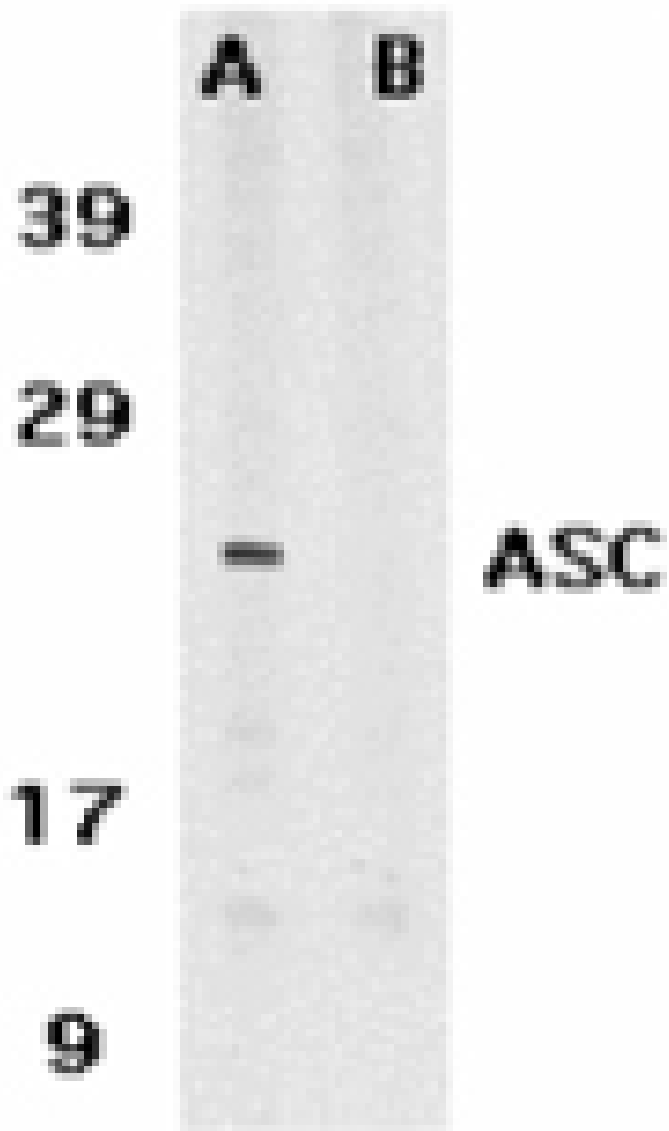
ADI-905-173-100	100µg
-----------------	-------

Manuals, SDS & CofA

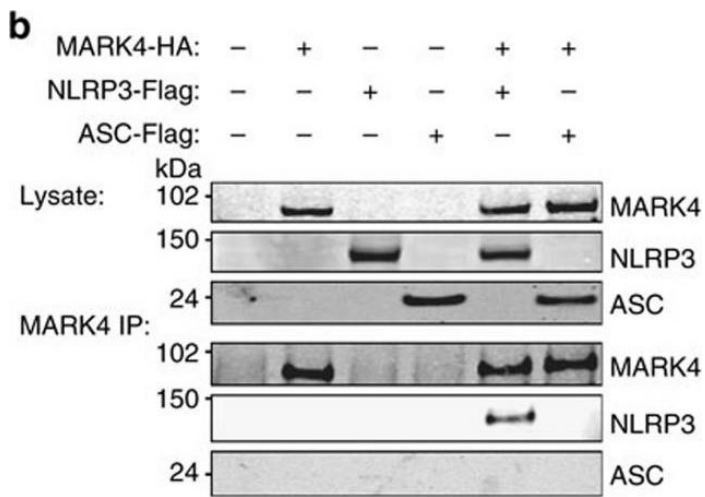
[View Online »](#)



Immunocytochemistry analysis of ASC in HL60 cells with ASC antibody at 5µg/ml.



Western blot analysis of ASC in HL60 whole cell lysate in the absence (A) or presence (B) of blocking peptide (2287P) with ASC antibody at 1µg/ml.



MARK4 interacts with NLRP3 in a microtubule-dependent manner. (a) Upon nigericin stimulation (3 μ M for 2 h), microtubule-disrupting drugs colchicine and nocodazole reduced the interaction between Mark4 and Nlrp3, shown by PLA signals in WT BMDM cells. Mark4 KO and Nlrp3 KO BMDM cells were employed as controls. (b) MARK4 was associated with NLRP3 in co-immunoprecipitation assay. Whole cell lysates were analysed as indication of transfection. Western blots are representative of 3 independent experiments. (c) Upon nigericin stimulation, PLA signal of NlrpP3 and Mark4, or Asc and Mark4 in BMDM cells derived from WT or Nlrp3 KO. Secondary only was employed as control in this PLA assay. (d) PLA signal of Nlrp3 and Asc in BMDM cells derived from WT or Nlrp3 KO or Asc KO before or after nigericin stimulation (3 μ M for 2 h). Mean \pm s.e.m. for all the cells taken from 5 to 8 different views at \times 40 magnification for each group (a,b,d). Comparisons of the two different groups were analysed by unpaired t-test. NS was considered as not statistically significant. *P<0.05, ***P<0.001 and ****P<0.0001 were considered as statistically significant (a,c,d). Results are representatives of three independent experiments. Scale bar, 10 μ m.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: MARK4 regulates NLRP3 positioning and inflammasome activation through a microtubule-dependent mechanism. *Nat Commun* (2017)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name

Apoptosis-associated speck-like protein containing a CARD, Caspase recruitment domain-containing protein 5, CARD-5, PYD and CARD domain-containing protein, PYCARD, Target of methylation-induced silencing, TMS1

Application

ICC, WB

Application Notes

Detects a band of ~25kDa by Western blot.

Formulation

Liquid. In PBS containing 0.02% sodium azide.

Host

Rabbit

Immunogen

Synthetic peptide corresponding to aa 182-195 of human ASC .

Purity Detail

Peptide affinity purified.

Recommendation Dilutions/Conditions

Western Blot (1µg/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

Q9ULZ3

Worry-free Guarantee

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

Last modified: May 29, 2024



Myo 212, Scd 124,
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-
fr@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)