

Caspase-11 (mouse) monoclonal antibody (8A5)

Caspase-11 not only activates caspase-1, that is required for the maturation of proinflammatory cytokines such as interleukin-1 (IL-1) and interleukin-18 (IL-18), but also activates caspase-3, leading to cellular apoptosis under pathological conditions. In most cells, caspase-11 is only expressed upon induction with pro-inflammatory stimuli. Caspase-11 is an essential mediator of endotoxic shock and caspase-11-deficient mice are resistant to endotoxic shock. LPS-induced caspase-11 regulates lymphocyte apoptosis by activating both caspase-3 and caspase-7. Human caspase-4, thought to be the homolog of mouse caspase-11, may be an effective therapeutic target for treatment of septic shock.

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

Ordering Information

[Order Online »](#)

ALX-804-494-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

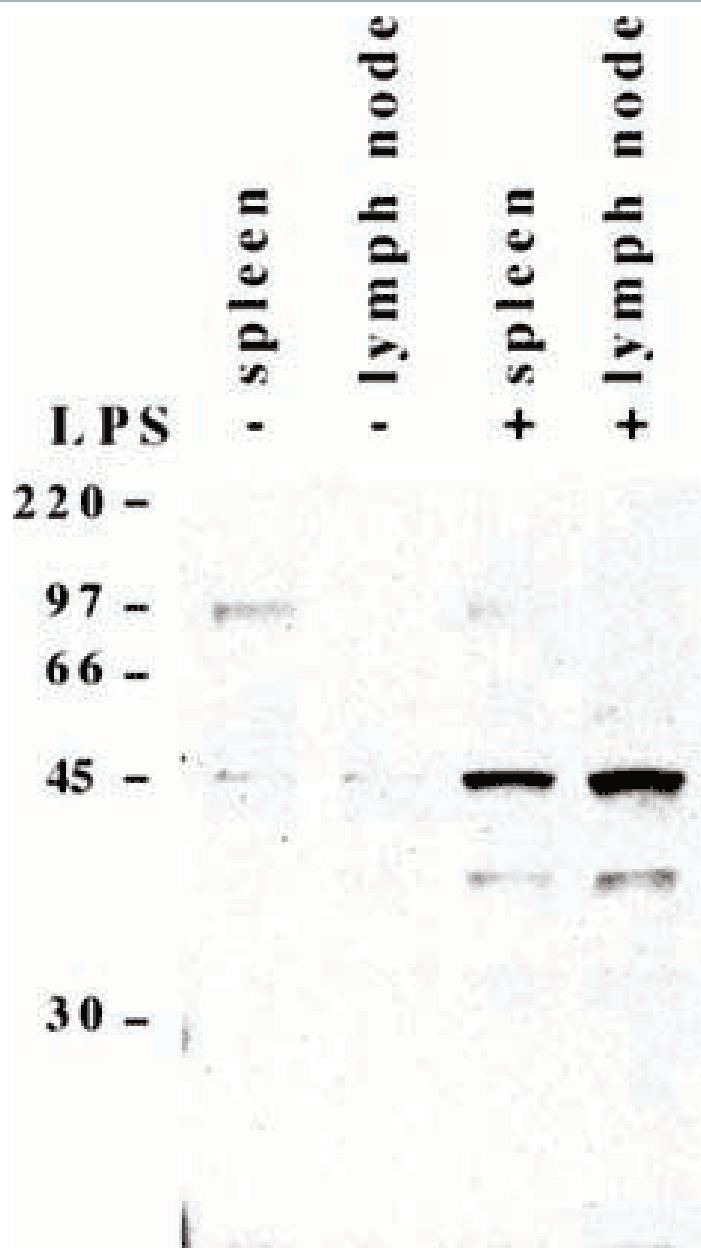


Figure: MAb 8A5 detects endogenous caspase-11 in mouse spleen and lymph node as two bands of 43 and 38 kDa. After 16 hrs co-culture with LPS 20 mg/mL elevated levels of caspase-11 are detected in mouse spleen and lymph node.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Flow Cytometry, WB

Application Notes Detects two bands of ~38kDa and ~43kDa by Western blot.

Clone 8A5

Formulation Liquid. In PBS containing 0.02% sodium azide.

Host Rat

Immunogen p20 fragment of mouse caspase-11.

Isotype IgG1

Purity Detail Protein G-affinity purified.

Species Reactivity Mouse, Rat

Specificity The epitope is in the p20 fragment.

UniProt ID P70343

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)