

# 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](#).

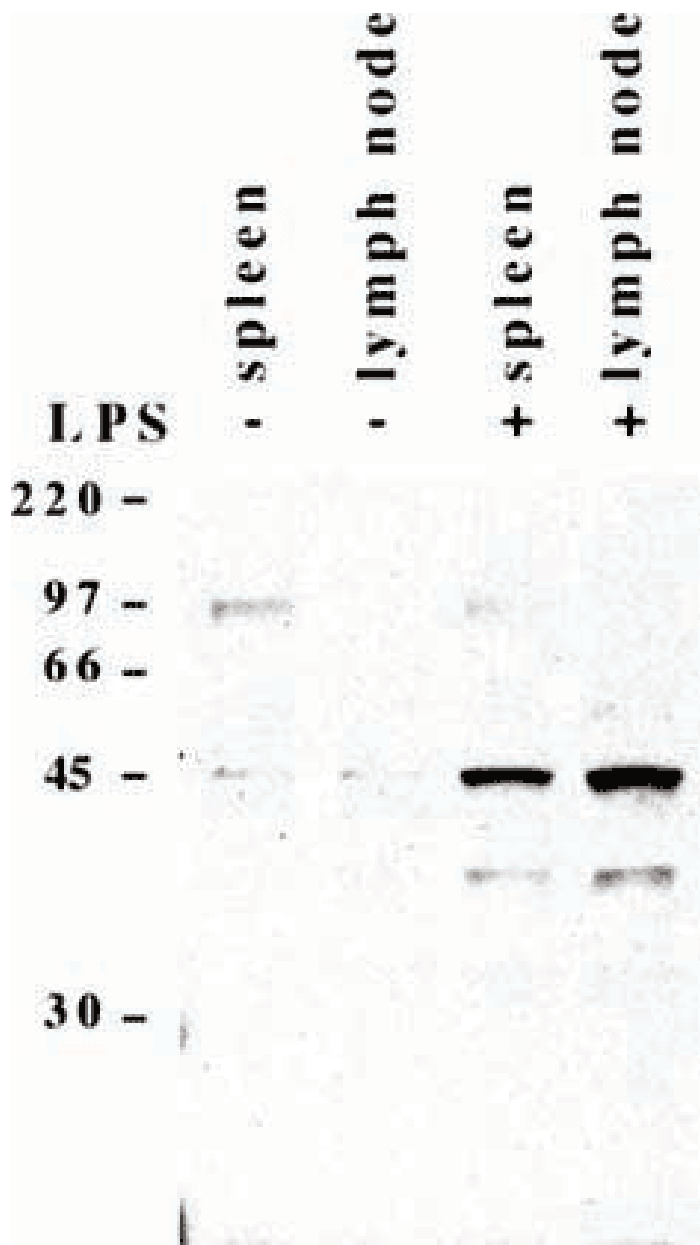
## Ordering Information

[Order Online »](#)

ALX-804-4940	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 <a href="#">Worry-Free Guarantee</a>

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-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)