

# FasL (mouse) monoclonal antibody (A11)

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

Citations: 11

[View Online »](#)

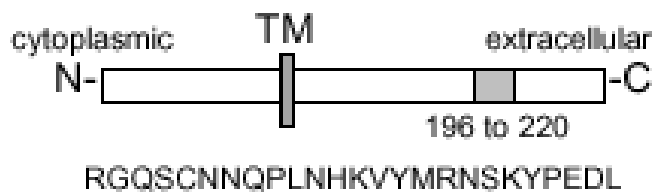
## Ordering Information

[Order Online »](#)

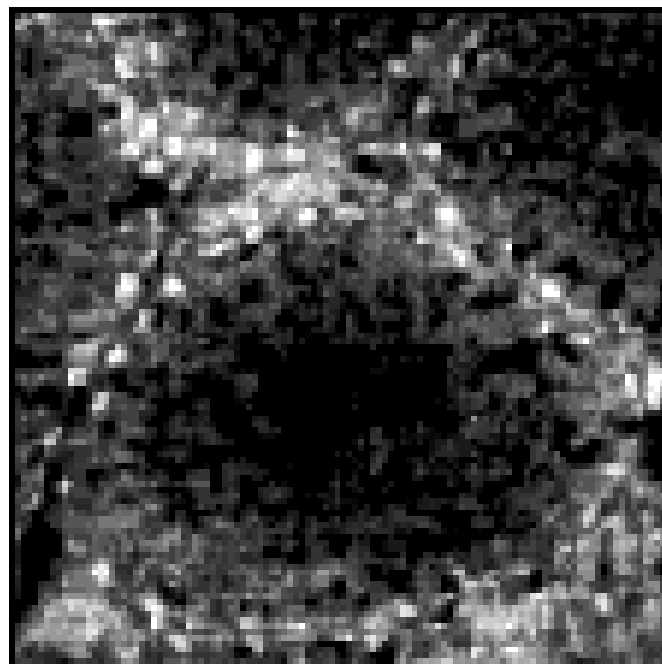
ALX-804-009-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Schematic structure of mouse FasL (peptide, aa. 196-220)



**Figure:** Immunohistochemical localization of FasL in adult mouse testis using the A11 antibody.

**Method:** Cryostat sections (6-8µm) were air dried, fixed in 4% paraformaldehyde in PBS and used immediately. Sections were incubated for 1 hr at room temperature (RT) with 20 µg/ml of A11 antibody in PBS, 0.2% BSA, 0.05% Tween 20. After rinsing in PBS, a FITC-conjugated anti-rat IgM was added for 30 min., slides washed again in PBS and visualized using a fluorescence microscope.

**Paraffin sections:** Dewax in xylene (2×10 min), EtOH 100% (2×10 min), block endogeneous peroxidase (0.6% H<sub>2</sub>O<sub>2</sub> in Methanol 10 min), rehydrate through graded EtOH (100%, 95%, 70%), rinse in H<sub>2</sub>O, unmask epitopes by heating slides in citrate buffer (0.01M, pH 6.0) 2×5 min in microwave oven (650-750W), cool down to RT 20 min, wash 1x in PBS, incubate with A11 20 µg/ml antibody, wash in PBS, incubate in anti-rat IgM HRP ([ALX-211-055](#)), wash in PBS, develop in appropriate substrate solution (DAB or AEC).", FasL (mouse) monoclonal antibody (A11) Immunohistochemistry, FasL (mouse) monoclonal antibody (A11) Immunohistochemistry, yes ALX-804-010, 804-

010.gif, <http://static.enzolifesciences.com/fileadmin/files/image/804-010.gif>, Schematic structure of mouse FasL (peptide

## Handling & Storage

Use/Stability	Stable for at least 6 months after receipt when stored at +4°C undiluted.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	+4°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	CD178, CD95L, APO-1L, TNFSF 6
Application	Flow Cytometry, IHC (FS), IHC (PS)
Application Notes	For Western blot, we recommend MAb to FasL (mouse) (H11) (Prod. No. ALX-804-010).
Clone	A11
Formulation	Liquid. Purified antibody in PBS containing 0.02% sodium azide.
Host	Rat
Immunogen	Synthetic peptide corresponding to aa 196-220 of mouse FasL (CD95L; APO-1L; CD178).
Isotype	IgM
Source	Purified from hybridoma tissue culture supernatant.
Species Reactivity	Mouse
UniProt ID	P41047
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a> .



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-  
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)