## FasL (mouse) monoclonal antibody (A11)

This antibody is covered by our Worry-Free Guarantee.

Citations: 11

View Online »

**Ordering Information** 

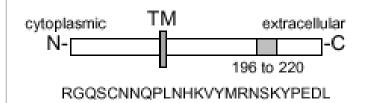
Order Online »

ALX-804-009-C100

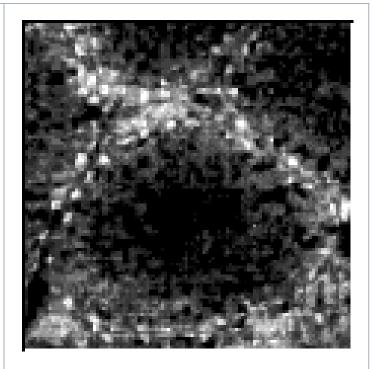
100µg

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/8 010.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 Worry-Free Guarantee

•

Last modified: May 29, 2024



info-

eu@enzolifesciences.com