

FasL (mouse) monoclonal antibody (H11)

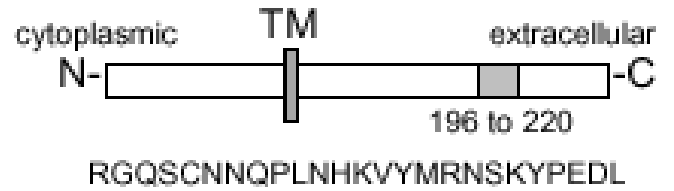
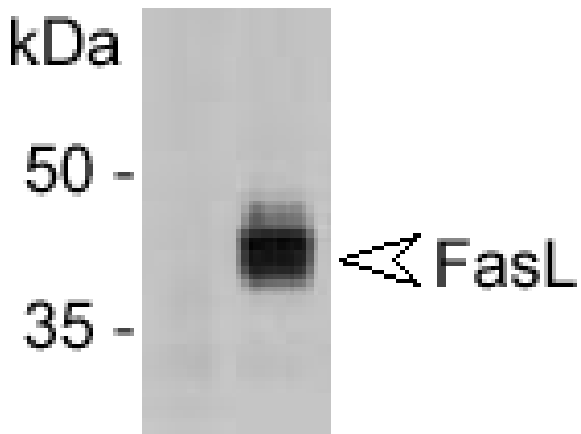
This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 4

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



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

Figure: Detection of FasL in 293T cells transfected with a FasL expression plasmid (right panel). Mock-transfected cells (left panel).

Method: Cells extracts from (2×10^6) cells transfected with a FasL-expression plasmid were resolved by SDS-PAGE under reducing conditions, transferred to nitrocellulose and probed with the monoclonal anti-FasL H11 antibody at $2\mu\text{g/ml}$. Proteins were visualized using a peroxidase-conjugated antibody to rat IgG ([ALX-211-052](#)) and a chemiluminescence detection system. Arrow indicates FasL. """, FasL (mouse) monoclonal antibody (H11) WB, FasL (mouse) monoclonal antibody (H11) Western blot, yes

ALX-804-015, 804-015_b.gif, http://static.enzolifesciences.com/fileadmin/files/image/804-015_b.gif, **Figure:** Immunolocalization of ryanodine receptor 1 (RyR1) in rat brain hippocampus using Monoclonal Antibody to Ryanodine Receptor (C3-33) (Prod. No. [ALX-804-015](#)). """"""

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, ICC, WB
Application Notes	For immunohistochemistry, we recommend MAb to FasL (A11) (Prod. No. ALX-804-009).
Clone	H11
Formulation	Liquid. 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	IgG2a
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](#)



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com

Last modified: May 29, 2024