FasL (mouse) monoclonal antibody (H11)

This antibody is covered by our Worry-Free Guarantee.

Citations: 4

View Online »

Ordering Information

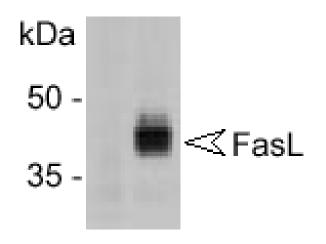
Order Online »

ALX-804-010-C100

Purified - 100µg

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 x 10⁶) 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 2μg/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.

Avoid freeze/thaw cycles. Handling

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).

lgG2a Isotype

Source Purified from hybridoma tissue culture supernatant.

Species Reactivity Mouse

UniProt ID P41047

Worry-free Guarantee

Last modified: May 29, 2024

This antibody is covered by our Worry-Free Guarantee



ENZO LIFE SCIENCES, Phone: 800.942.0430 info-

ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 usa@enzolifesciences.com

eu@enzolifesciences.com

European Sales Office

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

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

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

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