FasL monoclonal antibody (3C82)

This antibody is covered by our Worry-Free Guarantee.

Citations: 1

View Online »

Ordering Information

Order Online »

ALX-804-273-C100

100µg

Manuals, SDS & CofA

View Online »

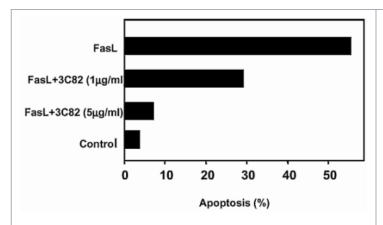
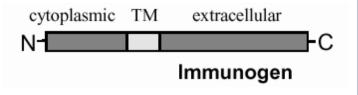


Figure: Inhibition of mouse FasL (supernatant of FasL-transfected Neuro2A cells) by 3C82.**Method:**Supernatants of Neuro2A cells containing murine FasL (1:2 dilution) were incubated with Fas ecpressing L929-APO-1 target cells. After 14h apoptosis was assessed by MTT staining.



Schematic structure of recombinant human FasL (extracellular domain).

Handling & Storage

Handling Do not freeze.

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name CD95L, APO-1L, CD178, TNFSF 6

Application Flow Cytometry, FUNC

Application Notes Functional Application: neutralizing

Clone 3C82

Formulation Liquid. In PBS, pH 7.2, containing 1% BSA. Contains no

preservatives.

Host Mouse

Immunogen Recombinant human FasL (APO-1L; CD95L;

CD178) (extracellular domain).

lgG1

Purity Detail Epitope-affinity purified.

Recommendation Dilutions/ConditionsSuggested starting dilution is >1:100Suggested

dilutions/conditions may not be available for all applications. Optimal conditions must be determined

individually for each application.

Species Reactivity Human, Mouse

UniProt ID P48023

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

•

Last modified: May 29, 2024



info-

eu@enzolifesciences.com