

FasL polyclonal antibody

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

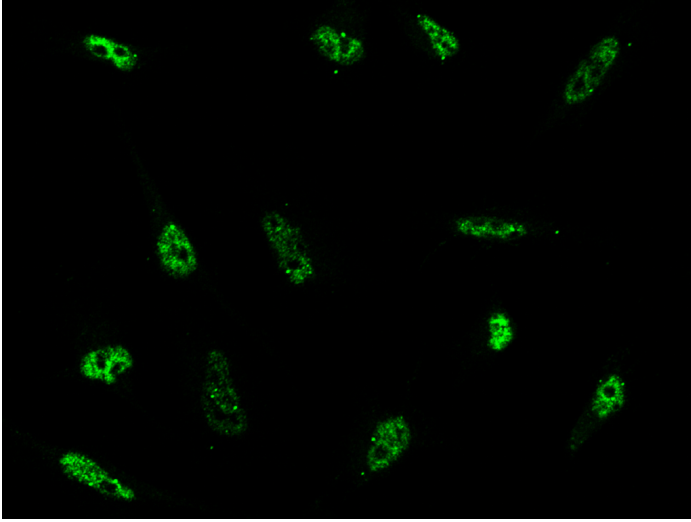
Ordering Information

[Order Online »](#)

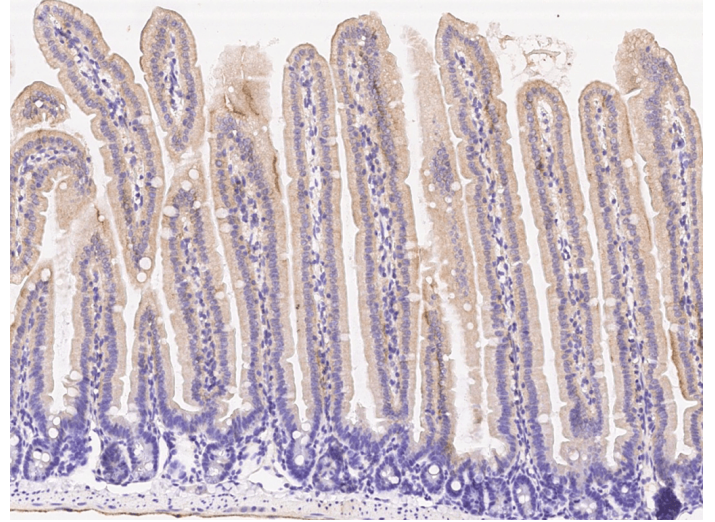
ENZ-ABS940-0100	100µl
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Immunofluorescence staining of FASLG in HeLa cells. Cells were fixed with 4% PFA, permeabilized with 0.3% Triton X-100 in PBS, blocked with 10% serum, and incubated with rabbit anti-human FASLG polyclonal antibody (1:1000) at 4 °C overnight. Then cells were stained with the Alexa Fluor®488-conjugated Goat Anti-rabbit IgG secondary antibody (green). Positive staining was localized to nucleus.



Immunochemical staining of mouse Faslg in mouse small intestine with rabbit polyclonal antibody (1:1000, formalin-fixed paraffin embedded sections).

Handling & Storage

Handling Protected from light. Avoid freeze-thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

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

Application ICC, IF, IHC (PS)

Formulation Liquid. 0.2 µm filtered solution in PBS. Preservative free.

Host Rabbit

Immunogen Synthetic peptide corresponding to the C-terminus of the mouse Fas Ligand / FASLG / CD95L.

Isotype IgG

Purity Detail Protein A and antigen affinity purified.

Recommendation Dilutions/Conditions Immunohistochemistry (paraffin sections) (1:500-1:2'000) Immunocytochemistry (1:300-1:10'000) Immunofluorescence (1:300-1:10'000) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Species Reactivity Mouse

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

Last modified: September 17, 2024



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