

CD81 polyclonal antibody

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

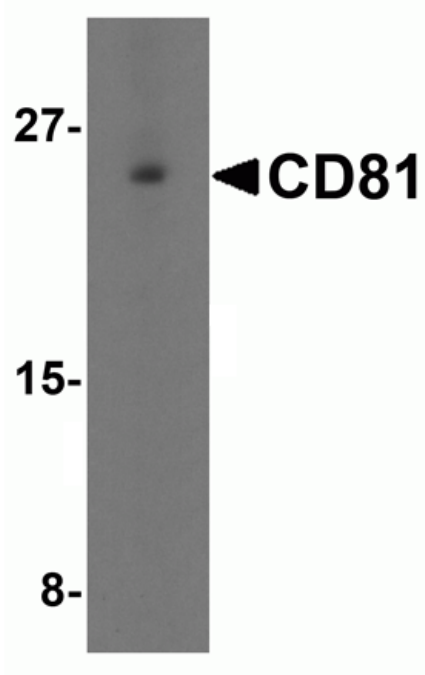
Ordering Information

[Order Online »](#)

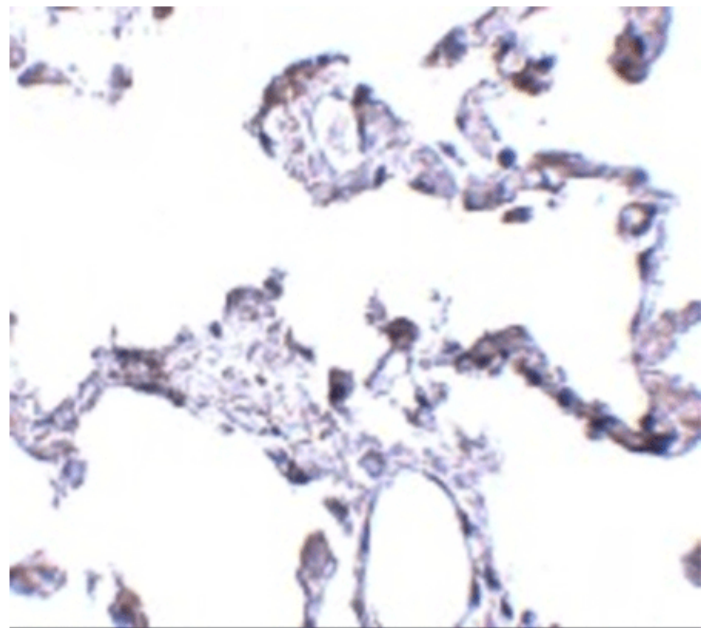
ENZ-ABS813-0100	100µg
-----------------	-------

Manuals, SDS & CofA

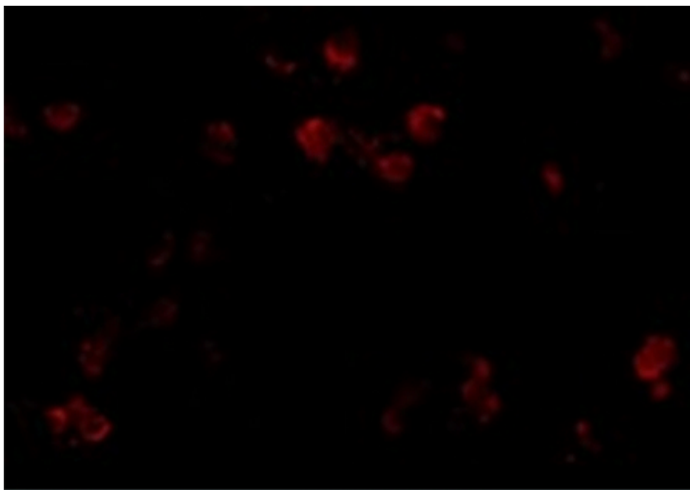
[View Online »](#)



Western blot analysis of CD81 in human lung tissue lysate with CD81 antibody at 1 $\mu\text{g/mL}$ in (A) the absence and (B) the presence of blocking peptide.



Immunohistochemistry of CD81 in human lung tissue with CD81 antibody at 5 $\mu\text{g/mL}$.



Immunofluorescence of CD81 in A549 cells with CD81 antibody at 20 $\mu\text{g/mL}$.

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 26 kDa cell surface protein TAPA-1, Tetraspanin-28, Tspan-28

Application ELISA, IF, IHC, WB

Formulation Liquid. In PBS containing 0.02% sodium azide

Host Rabbit

Immunogen The immunogen is located within amino acids 30 – 80 of CD81.

Isotype IgG

Purity Detail Affinity purified.

Species Reactivity Human, Mouse, Rat

UniProt ID P60033

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

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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