

DLL4 polyclonal antibody

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

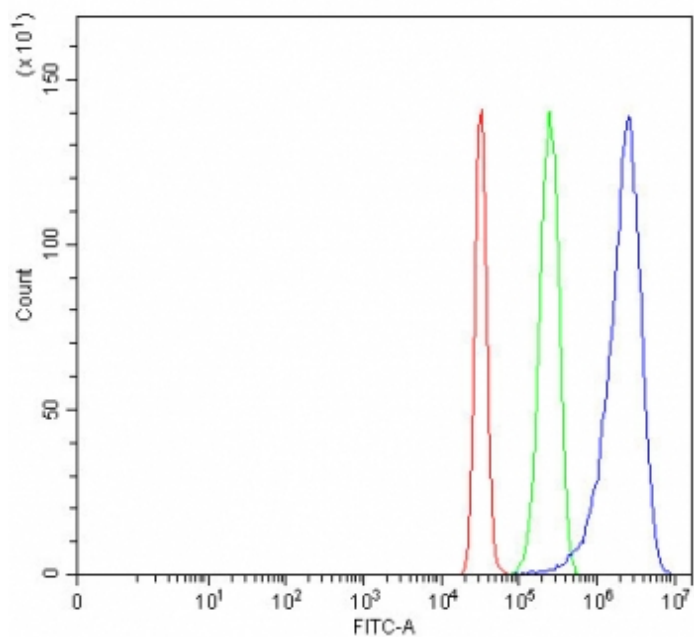
Ordering Information

[Order Online »](#)

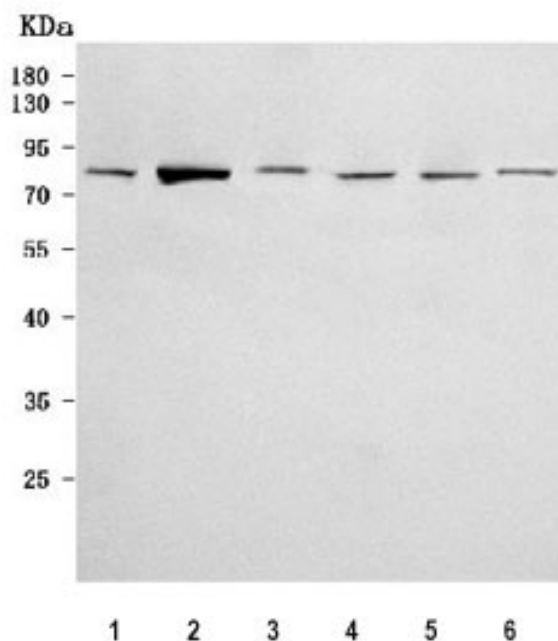
ENZ-ABS887-0100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Flow cytometry testing of fixed and permeabilized human HepG2 cells with Delta-like protein 4 antibody at $1\mu\text{g}/\text{million cells}$ (blocked with goat sera); Red=cells alone, Green=isotype control, Blue= Delta-like protein 4 antibody.



Western blot testing of 1) human A431, 2) human 293T, 3) human Caco-2, 4) human PC-3, 5) rat brain and 6) mouse brain tissue lysate with Delta-like protein 4 antibody. Predicted molecular weight $\sim 75\text{ kDa}$ but may be observed at higher molecular weights due to glycosylation.

Handling & Storage

Use/Stability	After reconstitution, the Delta-like protein 4 antibody can be stored for up to one month at 4°C. For long-term, aliquot and store at -20°C.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Delta-like protein 4
Application	ELISA, Flow Cytometry, WB
Formulation	Lyophilized from 1X PBS containing 2% Trehalose
Host	Rabbit
Immunogen	E. coli-derived recombinant human DLL4 (aa 27-678).
Isotype	IgG
Purity Detail	Antigen affinity purified.
Recommendation Dilutions/Conditions	ELISA (direct) (0.1-0.5µg/ml)Flow cytometry (1-3µg/million cells)Western blot (0.5-1µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse, Rat
UniProt ID	Q9NR61
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee .



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-
eu@enzolifesciences.com](mailto:info-eu@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)