## DLL4 polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

**Ordering Information** 

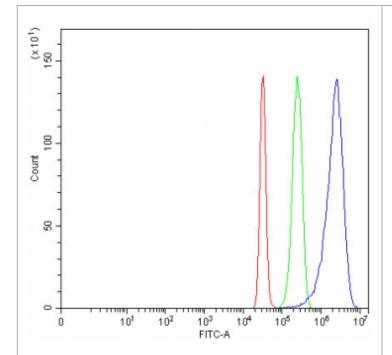
Order Online »

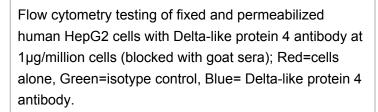
ENZ-ABS887-0100

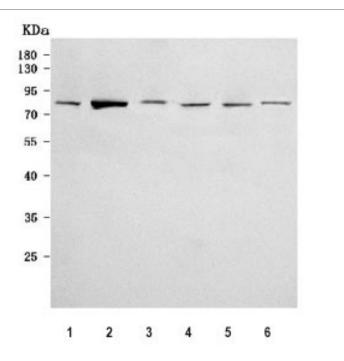
100µg

Manuals, SDS & CofA

**View Online** »







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 ~75 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

.

Last modified: July 12, 2024



info-

eu@enzolifesciences.com