

# DLL3 polyclonal antibody

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

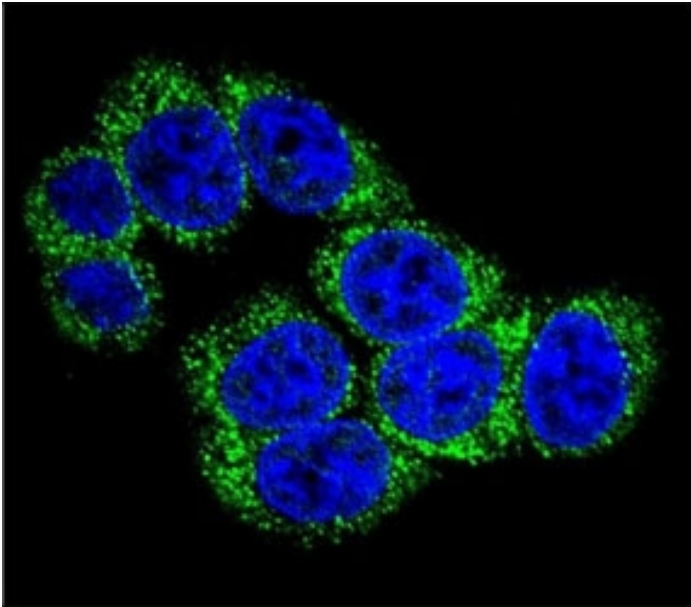
## Ordering Information

[Order Online »](#)

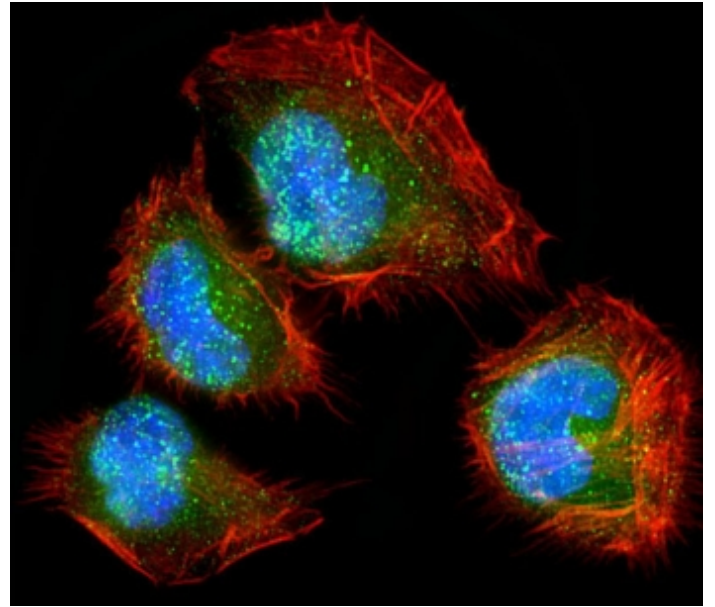
ENZ-ABS888-0200	0.2ml
-----------------	-------

## Manuals, SDS & CofA

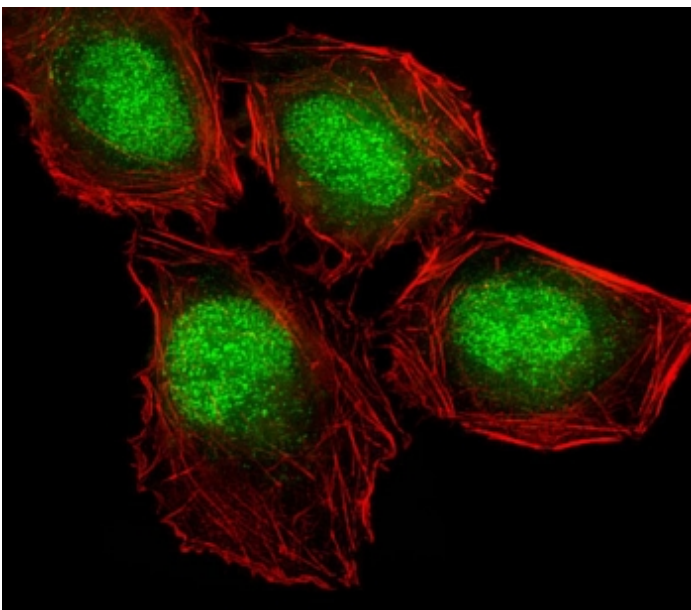
[View Online »](#)



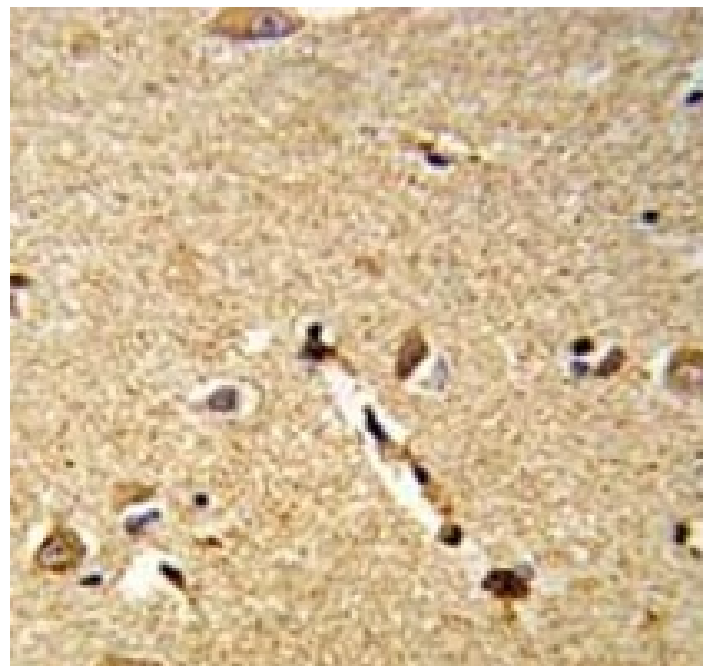
Confocal immunofluorescent analysis of DLL3 antibody with 293 cells followed by Alexa Fluor 488-conjugated goat anti-rabbit IgG (green). DAPI was used as a nuclear counterstain (blue).



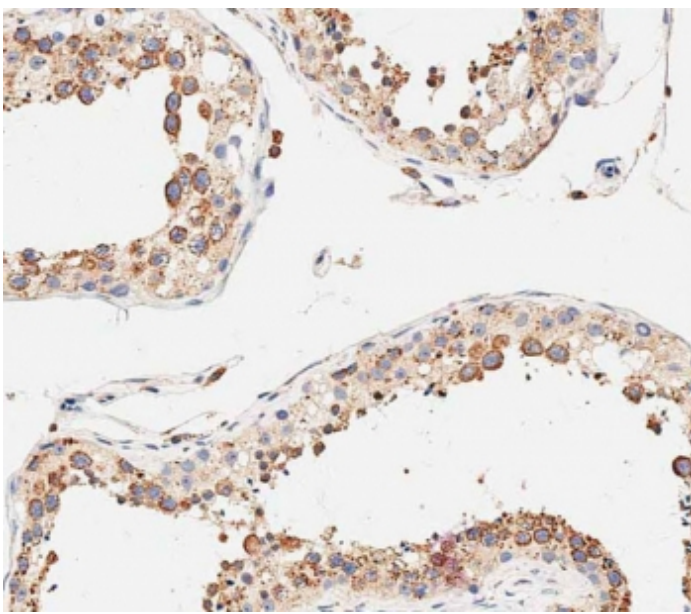
Immunofluorescent staining of human U-2 OS cells with DLL3 antibody (green), Phalloidin (red) and DAPI nuclear stain (blue).



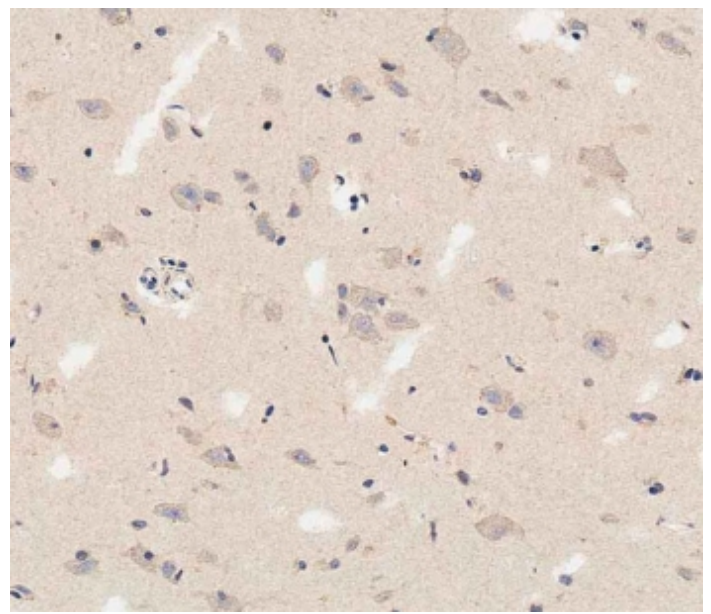
Immunofluorescent staining of human U-2 OS cells with DLL3 antibody (green) and Phalloidin (red).



IHC analysis of FFPE human brain tissue stained with DLL3 antibody



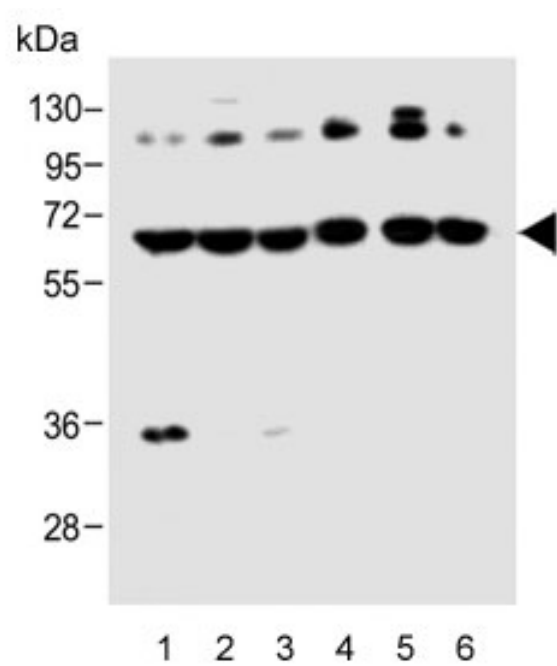
IHC staining of FFPE human testis with DLL3 antibody. HIER: boil tissue sections in pH 9 EDTA for 20 min and allow to cool before testing.



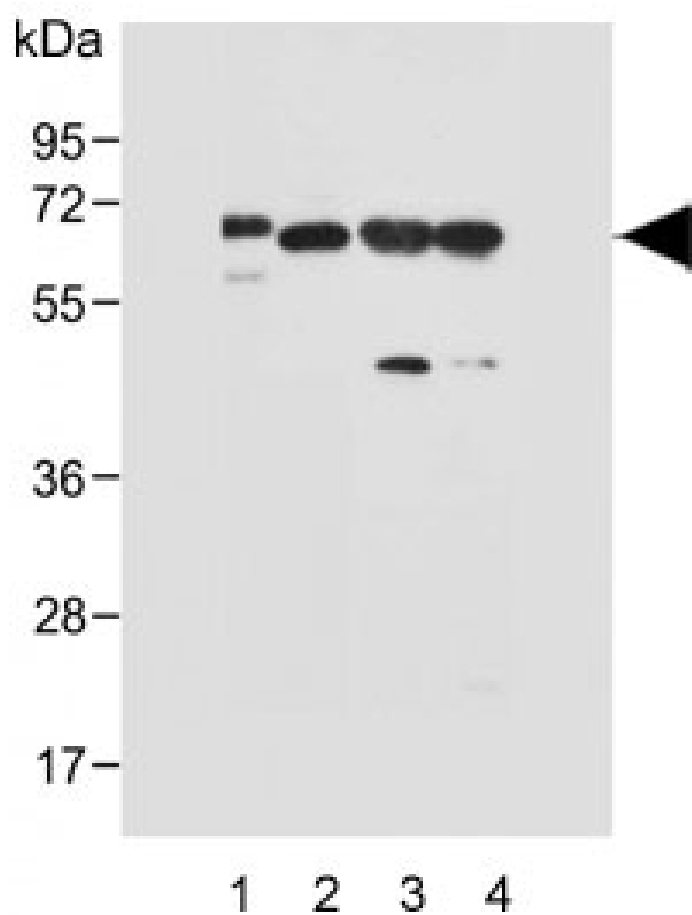
IHC staining of FFPE human brain with DLL3 antibody. HIER: boil tissue sections in pH 9 EDTA for 20 min and allow to cool before testing.



Western blot analysis of DLL3 antibody and A375 lysate. Predicted molecular weight: ~65 kDa.



Western blot testing of human 1) A375, 2) U-251 MG, 3) NCI-H1299, 4) brain, 5) mouse brain and 6) rat brain lysate with DLL3 antibody. Predicted molecular weight: ~65 kDa.



Western blot testing of human 1) NCI-H1299, 2) brain, 3) mouse brain and 4) rat brain lysate with DLL3 antibody. Predicted molecular weight: ~65 kDa.

## Handling & Storage

Handling	Aliquot the DLL3 antibody and store frozen at -20°C or colder. 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 3
Application	ELISA, IF, IHC, WB
Formulation	Liquid. In 1X PBS, pH 7.4, containing 0.09% sodium azide.
Host	Rabbit
Immunogen	Human DLL3 (aa 519-548).
Isotype	Ig
Purity Detail	Antigen affinity purified.
Recommendation Dilutions/Conditions	Immunofluorescence (1:10-1:50)Immunohistochemistry (paraffin sections) (1:50-1:100)Western blot (1:1,000)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	Q9NYJ7
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>