

Polysialic acid Recombinant monoclonal antibody (735) (Rabbit IgGκ)

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

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-ABS559-0100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Stable at 4°C for up to 3 months. For longer storage, aliquot and store at -20°C.
Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	PSA, N-CAM, Neural cell adhesion molecule, polySA
Application	Flow Cytometry, IHC, IP
Clone	735
Formulation	Liquid. In PBS containing 0.02% Proclin 300.
Host	Rabbit
Immunogen	Homopolymers of α 2,8-linked sialic acid of <i>E. coli</i> K1.
Isotype	IgGk
Purity Detail	Protein A affinity purified.
Species Reactivity	Human, Mouse, Rat
Specificity	Recognizes polysialic acid (PSA), a carbohydrate polymer attached to the neural cell adhesion molecule (NCAM), which shares molecular epitopes with homopolymers of α 2,8-linked sialic acid of <i>E. coli</i> K1.

Technical Info / Product Notes This chimeric rabbit antibody was made using the variable domain sequences of the original murine IgG2a format, for improved compatibility with existing reagents, assays and techniques.



ENZO LIFE SCIENCES
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
& Luxembourg
(ELS) AG
Phone: +41 61 926 8989
info-be@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@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

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee.