Polysialic acid Recombinant monoclonal antibody (735) (Mouse IgG2aκ)

This antibody is covered by our Worry-Free Guarantee.

Citations: 1

View Online »

Ordering Information

Order Online »

ENZ-ABS560-0200

200µg

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 Mouse

Immunogen Homopolymers of α2,8-linked sialic acid of E. coli K1.

Isotype IgG2ak

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

 $\alpha 2,8\text{-linked}$ sialic acid of \textit{E. coli} K1.

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

•

Last modified: May 29, 2024



info-

eu@enzolifesciences.com