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

lgGk

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

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

Technical Info / Product NotesThis 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.

Last modified: May 29, 2024



info-

eu@enzolifesciences.com