CD56 Recombinant monoclonal antibody (RM315)

CD56, also called NCAM1 and Neural cell adhesion molecule 1, is a cell adhesion molecule involved in neuron-neuron adhesion, neurite fasciculation, outgrowth of neurites, etc. Acts as a receptor for rabies virus.

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS673-0100

100µl

Manuals, SDS & CofA

View Online »

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Neural cell adhesion molecule 1, NCAM-1

Application IHC (PS), WB

Clone RM315

Formulation Liquid. In 1x PBS containing 50% glycerol, 1% BSA and

0.09% sodium azide

Host Rabbit

Immunogen Synthetic peptide corresponding to the C-terminus of

human CD56.

Isotype IgG

Purity Detail Protein A affinity purified.

Recommendation Dilutions/Conditions Immunohistochemistry (FFPE): 1:500-1:1,000Western

Blot: 1:1,000-1:2,000Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Source Purified from animal origin-free supernatant.

Species Reactivity Human

UniProt ID P13591

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

•

Last modified: May 29, 2024



info-

eu@enzolifesciences.com