CD31 Recombinant monoclonal antibody (RMCD31-1)

CD31 is found on the surface of platelets, monocytes, neutrophils, and some types of T-cells, and makes up a large portion of endothelial cell intercellular junctions. It is a member of the immunoglobulin superfamily and is likely involved in leukocyte migration, angiogenesis, and integrin activation. [RefSeq]

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS340-0100

100µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name PECAM-1, Platelet endothelial cell adhesion molecule

Application IF, IHC (PS)

Clone RMCD31-1

Formulation Liquid. In 1X PBS containing 0.1 mg/ml BSA (US sourced)

and 0.05% sodium azide.

Gene/Protein Identifier 5175 (Gene ID)

Host Rabbit

Immunogen Recombinant full length human CD31.

Isotype IgGk

Purity Detail Protein A affinity purified.

Recommendation Dilutions/Conditions Immunofluorescence (0.5-1 µg/mL) Immunohistochemistry

(formalin-fixed paraffin sections) (0.5-1 μ g/mL) for 30 minutes at RTSuggested dilutions/conditions may not be available for all applications.Optimal conditions must be

determined individually for each application.

Species Reactivity Human

UniProt ID P16284

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

•

Last modified: May 29, 2024



info-

eu@enzolifesciences.com