CD34 Recombinant monoclonal antibody (My10) (Rabbit IgGк)

Human CD34, an \sim 105 - 120 kDa type 1 transmembrane glycoprotein expressed on the surface of most human hematopoietic stem and progenitor cells (HSPCs) as well as on mesenchymal stem cells, embryonic fibroblasts, endothelial cells, neurons, and some tumor cell lines.

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS628-0200 20

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 Hematopoietic progenitor cell antigen CD35

Application Flow Cytometry, IHC, IP

Clone My10

Formulation Liquid. In PBS containing 0.02% Proclin 300.

Host Rabbit

Immunogen KG-1a Human myeloblastic leukemic cell line (immature

human myeloid cell line) from a patient with

nonlymphocytic leukemia.

Isotype IgGk

Purity Detail Protein A affinity purified.

Species Reactivity Human

Technical Info / Product NotesThis chimeric rabbit antibody was made using the variable

domain sequences of the original murine IgG1 format, for improved compatibility with existing reagents, assays and

techniques.

UniProt ID P28906

Last modified: May 29, 2024



info-

eu@enzolifesciences.com