CD20 Recombinant monoclonal antibody (RMC20-1)

CD20 is expressed on all stages of B cell development except the first and last; it is present from late pro-B cells through memory cells, but not on either early pro-B cells or plasma blasts and plasma cells. It has no known natural ligand and its function is to enable optimal B-cell immune response, specifically against T-independent antigens. It is suspected that it acts as a calcium channel in the cell membrane. It has been shown that CD20 plays a role in the microenvironmental interactions of B cells. [Wiki]

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS346-0100 10

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 B-lymphocyte antigen CD20

Application IHC (PS)

Clone RMC20-1

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

and 0.05% sodium azide.

Host Rabbit

Immunogen Full length human CD20.

Isotype IgGk

Purity Detail Protein A affinity purified.

Recommendation Dilutions/Conditions Immunohistochemistry (formalin-fixed paraffin sections)

(0.5-1 µg/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 P11836

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

.



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com