CD20 Recombinant monoclonal antibody (10F381) (Human IgG1ĸ)

CD20 is a hydrophobic transmembrane protein with a molecular weight of approximately 35 kDa located on pre-B and mature B lymphocytes, and most B-cell non-Hodgkin's lymphomas, but is not found on stem cells, pro-B cells, normal plasma cells or other normal tissues.

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS419-0200

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 B-lymphocyte antigen CD20, B-lymphocyte surface

antigen B1, Bp35 Leukocyte surface antigen Leu-16, Membrane-spanning 4-domains subfamily A member 1

Application Flow Cytometry

Clone 10F381 (Rituximab)

Formulation Liquid. In PBS containing 0.02% Proclin 300.

Host Human

Immunogen Human lymphoblastoid cell line SB.

lgG1κ

Purity Detail Protein A affinity purified.

Species Reactivity Human

UniProt ID P11836

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 infousa@enzolifesciences.com 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

Phone: +49 7621 5500 526 infode@enzolifesciences.com UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com