

CD20 Recombinant monoclonal antibody (10F381) (Rabbit IgGκ)

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-ABS418-0200	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	Rabbit
Immunogen	Human lymphoblastoid cell line SB.
Isotype	IgG _k
Purity Detail	Protein A affinity purified.
Species Reactivity	Human
Technical Info / Product Notes	This chimeric rabbit antibody was made using the variable domain sequences of the original Human IgG1 format, for improved compatibility with existing reagents, assays and techniques.
UniProt ID	P11836

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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