



PRODUCT DATA SHEET

CD20 Recombinant monoclonal antibody (rIGEL/773)

ENZ-ABS316

Product Number/Sizes

ENZ-ABS316-0100 100 µg

Product Specifications

ALTERNATIVE NAME: B-lymphocyte antigen CD20
CLONE: rIGEL/773
HOST: Mouse
ISOTYPE: IgG2aκ
IMMUNOGEN: Recombinant full length human CD20.
UNIPROT ID: P11836
SPECIES REACTIVITY: Human
APPLICATIONS: Flow Cytometry, IF, IHC (PS)
RECOMMENDED DILUTIONS/CONDITIONS: Flow Cytometry (0.5-1 µg/10⁶ cells in 0.1 mL)
Immunofluorescence (0.5-1 µg/mL)
Immunohistochemistry (formalin-fixed paraffin sections) (0.5-1 µg/mL) for 30 minutes at RT
Suggested dilutions/conditions may not be available for all applications.
Optimal conditions must be determined individually for each application.
APPLICATION NOTES: This mAb can be used for immunophenotyping of leukemia and malignant cells, B lymphocyte detection in peripheral blood and B cell localization in tissues.
PURITY DETAIL: Protein G affinity purified.
FORMULATION: Liquid. In 1X PBS containing 0.1 mg/ml BSA (US sourced) and 0.05% sodium azide.
SHIPPING: Shipped on Blue Ice
LONG TERM STORAGE: +4°C
SCIENTIFIC BACKGROUND: CD20 is expressed by pre B-cells and persists during all stages of B-cell maturation but is lost upon terminal differentiation into plasma cells. It reacts with the majority of B-cells present in peripheral blood and lymphoid tissues and their derived lymphomas. In lymphoid tissue, germinal center blasts and B-immunoblasts are particularly reactive. It is a reliable antibody for ascribing a B-cell phenotype in known lymphoid tissues. Rarely, CD20-positive T-cell lymphomas have been reported. Reactivity has also been noted with Reed-Sternberg cells in cases of Hodgkin s disease, particularly of lymphocyte predominant type.

Revised 20-Dec-18

For Research Use Only, Not for Human

GLOBAL HEADQUARTERS

Enzo Life Sciences, Inc.
10 Executive Blvd
Farmingdale, NY 11735
USA
T 1-800-942-0430
T 1-631-694-7070
F 1-610-941-9252
E info-usa@enzolifesciences.com
www.enzolifesciences.com

EUROPE/ASIA

Enzo Life Sciences (ELS) AG
Industriestrasse 17, Postfach
CH-4415 Lausen
Switzerland
T +41/061 926 89 89
F +41/061 926 89 79
E info-ch@enzolifesciences.com
www.enzolifesciences.com