



# PRODUCT DATA SHEET

## CD20 Recombinant monoclonal antibody (IGEL/1497R)

ENZ-ABS345

### Product Number/Sizes

ENZ-ABS345-0100 100 µg

### Product Details

ALTERNATIVE NAME:	B-lymphocyte antigen CD20
CLONE:	IGEL/1497R
HOST:	Rabbit
ISOTYPE:	IgGκ
IMMUNOGEN:	Recombinant human full-length CD20.
UNIPROT ID:	P11836
SPECIES REACTIVITY:	Human
APPLICATIONS:	Flow Cytometry, IF, IHC (PS)
RECOMMENDED DILUTIONS/CONDITIONS:	Flow Cytometry (0.5-1 µg/10 <sup>6</sup> cells) 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.
PURITY DETAIL:	Protein A affinity purified.
FORMULATION:	Liquid. In 1X PBS containing 0.1 mg/ml BSA (US sourced) and 0.05% sodium azide.
SHIPPING:	Blue Ice
LONG TERM STORAGE:	+4°C
SCIENTIFIC BACKGROUND:	CD20 is a non-Ig differentiation antigen of B-cells and its expression is restricted to normal and neoplastic B-cells, being absent from all other leukocytes and tissues. It is expressed by pre B-cells and persists during all stages of maturation but is lost upon terminal differentiation into plasma cells.
REGULATORY STATUS:	RUO - Research Use Only

Revised 19-Dec-23

#### 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](mailto:info-usa@enzolifesciences.com)  
[www.enzolifesciences.com](http://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](mailto:info-ch@enzolifesciences.com)  
[www.enzolifesciences.com](http://www.enzolifesciences.com)