

CD20 Recombinant monoclonal antibody (IGEL/1497R)

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.

This antibody is covered by our [Worry-Free Guarantee](#).

Ordering Information

[Order Online »](#)

| | |
|-----------------|-------|
| ENZ-ABS345-0100 | 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 | Flow Cytometry, IF, IHC (PS) |
| Clone | IGEL/1497R |
| Formulation | Liquid. In 1X PBS containing 0.1 mg/ml BSA (US sourced) and 0.05% sodium azide. |
| Host | Rabbit |
| Immunogen | Recombinant human full-length CD20. |
| Isotype | IgGκ |
| Purity Detail | Protein A affinity purified. |
| Recommendation Dilutions/Conditions | Flow Cytometry (0.5-1 µg/106 cells)Immunofluorescence (0.5-1 µg/mL)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](#)



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-](mailto:info-usa@enzolifesciences.com)
usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-](mailto:info-fr@enzolifesciences.com)
fr@enzolifesciences.com

France
Phone: +33 472 440 655
[info-](mailto:info-fr@enzolifesciences.com)
fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
[info-](mailto:info-de@enzolifesciences.com)
de@enzolifesciences.com

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

Last modified: May 29, 2024