

CD22 Recombinant monoclonal antibody (hL22) (Rabbit IgGκ)

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

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-cell receptor CD22, B-lymphocyte cell adhesion molecule, BL-CAM, Sialic acid-binding Ig-like lectin 2, Siglec-2, T-cell surface antigen Leu-14, Epratuzumab
Application	Flow Cytometry, FUNC, IP
Clone	hL22
Formulation	Liquid. In PBS containing 0.02% Proclin 300.
Host	Rabbit
Immunogen	This antibody was prepared by the humanization of LL2 (EPB-2), a murine anti-CD22 IgG2a raised against Raji Burkitt lymphoma cells. Murine sequences comprise 5–10% of the molecule, with the remainder being human framework sequences, which greatly reduces the potential for immunogenicity.
Isotype	IgGk
Purity Detail	Protein A affinity purified.
Species Reactivity	Cynomolgus, Human, Rhesus monkey
Specificity	Recognizes the 3rd Ig-like domain of human CD22 (epitope B), a cell surface glycoprotein present on mature B-cells and on many types of malignant B-cells.

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

P20273

Worry-free Guarantee

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

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