CD22 Recombinant monoclonal antibody (hL22) (Rabbit IgGк)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS620-0200

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-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



info-

eu@enzolifesciences.com