CD22 Recombinant monoclonal antibody (hL22) (Human IgG1κ)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS619-0200

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 Human

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.

lgG1κ

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.

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



info-

eu@enzolifesciences.com