CD23 (human) monoclonal antibody (9p25)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS248-0100

100µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Short Term Storage -20°C

Long Term Storage -80°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Fc-ε-RII, BLAST-2

Application Flow Cytometry, FUNC

Application Notes Functional Application: Study of IgE binding to B cells

Clone 9p25

Formulation Liquid. In TRIS-NaCl, pH 8.

Host Mouse

Immunogen EBV transformed lymphoblastoid cell line.

lsotype lgG1

Purity Detail Purified using QMA Hyper D ion exchange

chromatography

Species Reactivity Human

Specificity Recognizes human B lymphocytes.

UniProt ID P06734

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg
Phone: +32 3 466 0420
infobe@enzolifesciences.com

Phone: +33 472 440 655 infofr@enzolifesciences.com Germany Phone: +49 7621 5500 526 infode@enzolifesciences.com UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com