CD19 Recombinant monoclonal antibody (4G7) (Mouse IgG1k)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS413-0200

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-lymphocyte antigen CD19, B-lymphocyte surface

antigen B4, Differentiation antigen CD19 T-cell surface

antigen Leu-12

Application Flow Cytometry, IF

Clone 4G7

Formulation Liquid. In PBS containing 0.02% Proclin 300.

Host Mouse

Immunogen Human CCL (chronic lymphocytic leukemia) cells.

IgG1ĸ

Purity Detail Protein A affinity purified.

Species Reactivity Human

UniProt ID P15391

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 infousa@enzolifesciences.com

ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
infoeu@enzolifesciences.com

European Sales Office

Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

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