CD30 Recombinant monoclonal antibody (Ki-1/1505R)

CD30 is the receptor for TNFSF8/CD30L. May play a role in the regulation of cellular growth and transformation of activated lymphoblasts. Regulates gene expression through activation of NF-kappa-B. [UniProt]

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS354-0100

100µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Ki-1, TNFRSF 8

Application Flow Cytometry, IF, IHC (PS)

Clone Ki-1/1505R

Formulation Liquid. In 1X PBS containing 0.1 mg/ml BSA (US sourced)

and 0.05% sodium azide.

Host Rabbit

Immunogen Partial human CD30.

Isotype IgGk

Purity Detail Protein A affinity purified.

Recommendation Dilutions/Conditions Flow Cytometry (0.5-1 µg/106 cells in 0.1

mL)Immunofluorescence (1-2

μg/mL)Immunohistochemistry (formalin-fixed paraffin sections) (0.5-1 μg/mL for 30 min at RT)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined

individually for each application.

Species Reactivity Human

UniProt ID P28908

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee



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

France
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

Last modified: May 29, 2024