CD30 Recombinant monoclonal antibody (Ki-3) (Mouse IgG2bк)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS432-0200

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 TNFRSF8, Tumor necrosis factor receptor superfamily

member 8, CD30L receptor, Ki-1 antigen, Lymphocyte

activation antigen CD30

Application ELISA, Flow Cytometry, IF, IHC

Application Notes Functional Application: Activation

Clone Ki-3

Formulation Liquid. In PBS containing 0.02% Proclin 300.

Host Mouse

Immunogen Human CD30.

lgG2bк

Purity Detail Protein A affinity purified.

Species Reactivity Human

UniProt ID P28908

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

.

Last modified: May 29, 2024



info-

eu@enzolifesciences.com