CD55 Recombinant monoclonal antibody (LU30) (Mouse IgG1λ)

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ENZ-ABS629-0200

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 Complement decay-accelerating factor, DAF

Application ELISA, IHC, WB

Clone LU30

Formulation Liquid. In PBS containing 0.02% Proclin 300.

Host Mouse

Immunogen Human CD55.

lgG1λ

Purity Detail Protein A affinity purified.

Species Reactivity Human

Specificity The antibody binds to CD55 with a Kd of 13 nM.

Technical Info / Product NotesThis chimeric mouse antibody was made using the

variable domain sequences of the original Human format, for improved compatibility with existing reagents, assays

and techniques.

UniProt ID P08174

Last modified: May 29, 2024



info-

eu@enzolifesciences.com