

# CD29 Recombinant monoclonal antibody (K20) (Mouse IgG2ak)

This antibody is covered by our [Worry-Free Guarantee](#).

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	Integrin $\beta$ -1
Application	Flow Cytometry, IHC (FS)
Clone	K20
Formulation	Liquid. In PBS containing 0.02% Proclin 300.
Host	Mouse
Immunogen	Human T-cell lymphoma cells injected into Biozzi high-responder mice.
Isotype	IgG2ak
Purity Detail	Protein A affinity purified.
Species Reactivity	Human
UniProt ID	P05556

# Worry-free Guarantee

This antibody is covered by our [Worry-Free Guarantee](#)