LRRC32 monoclonal antibody (Plato-1) (FITC conjugate)

LRRC32 (leucine rich repeat containing 32; also known as GARP or Garpin; Glycoprotein A repetitions predominant) is a glycoprotein expressed on the cell surface of megakaryocytes, platelets and activated regulatory T (Treg) cells.

This antibody is covered by our Worry-Free Guarantee.

Citations: 6

View Online »

Ordering Information

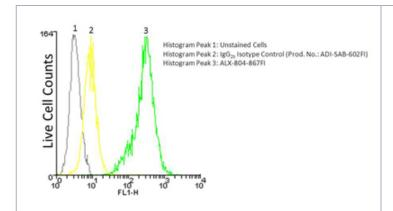
Order Online »

ALX-804-867FI-0100

100µg

Manuals, SDS & CofA

View Online »



Flow cytometry analysis, intracellular staining, of 10^6 Jurkat cells using LRRC32, mAb (Plato-1) (FITC conjugate) at a concentration of $50\mu g/ml$.

Handling & Storage

Use/Stability Stable for at least 1 year after receipt when stored at -20°C.

Handling Keep conjugated formats at +4°C. Do not freeze. Protect from light.

Short Term Storage +4°C

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Leucine-rich repeat-containing protein 32, GARP, Garpin,

Glycoprotein A repetitions predominant

Application Flow Cytometry

Application NotesDetects a band of ~100kDa by Western blot. Calculated

molecular weight of the full-length protein is ~72kDa.

Clone Plato-1

Formulation Liquid. In PBS containing 0.09% sodium azide.

Host Mouse

Immunogen Recombinant human LRRC32 (aa 20-627).

Isotype IgG2b

Purity Detail Protein G-affinity purified.

Recommendation Dilutions/Conditions Flow Cytometry (1:20)Suggested dilutions/conditions may

not be available for all applications. Optimal conditions must be determined individually for each application.

Source Purified from concentrated hybridoma tissue culture

supernatant.

Species Reactivity

Human, Mouse

UniProt ID

Q14392

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



info-

ENZO LIFE SCIENCES, European Sales Office ENZO LIFE SCIENCES Phone: 800.942.0430 (ELS) AG Phone: +41 61 926 8989 usa@enzolifesciences.com eu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

Phone: +33 472 440 655 fr@enzolifesciences.com

Phone: +49 7621 5500 526 de@enzolifesciences.com

UK & Ireland Phone (UK customers): 0845 601 1488 Phone: +44 1392 825900 info-uk@enzolifesciences.com