

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.

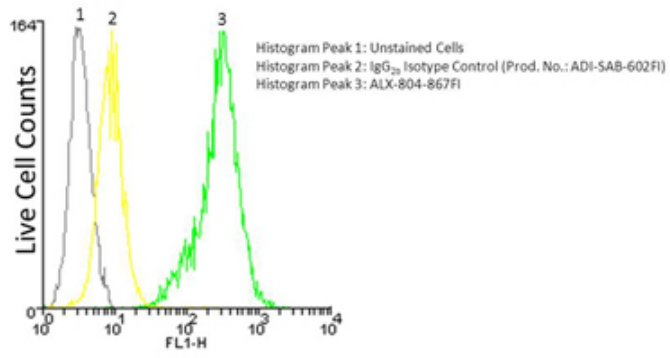
This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 6

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Flow cytometry analysis, intracellular staining, of 10⁶ Jurkat cells using LRRC32, mAb (Plato-1) (FITC conjugate) at a concentration of 50µg/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 Notes	Detects 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



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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