

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 »](#)

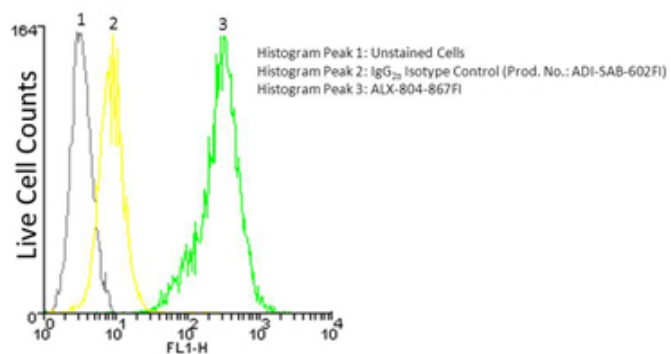
Ordering Information

[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 μ 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 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](#).



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)