

# LRRC32 monoclonal antibody (Plato-1) (PE 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: 3

[View Online »](#)

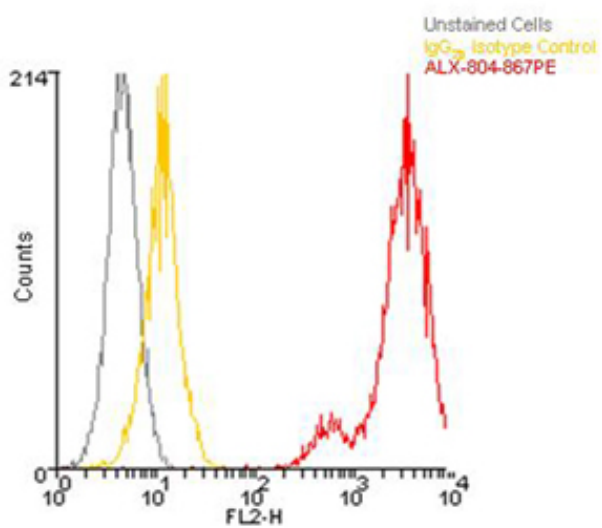
Ordering Information

[Order Online »](#)

ALX-804-867PE-0100	100µg
--------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Flow cytometry analysis, intracellular stained, of  $10^6$  Jurkat cells using LRRC32, mAb (Plato-1) (PE conjugate) at a concentration of  $10\mu\text{g/ml}$ .

## 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. Native 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 (1:100)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)