

LRRC32 (human):Fc (human), (recombinant)

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 (T_{reg}) cells.

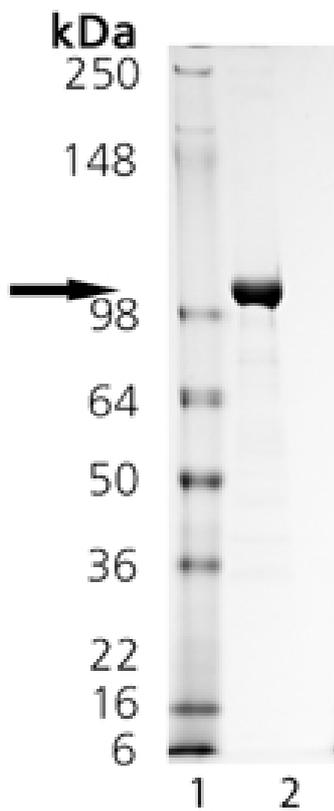
Ordering Information

[Order Online »](#)

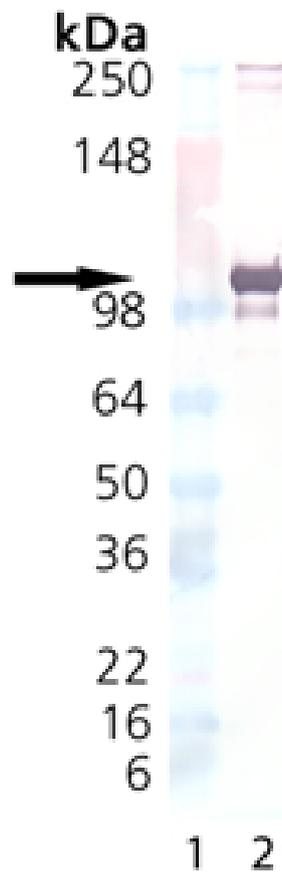
ALX-522-117-C050	50µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE analysis: Lane 1: MW Marker, Lane 2: 2 μ g LRRC32 (human):Fc (human), (recombinant).



Western Blot analysis: Lane 1: MW Marker, Lane 2: 100ng LRRC32 (human):Fc (human), (recombinant) probed with LRRC32 mAb (Prod. no. ALX-804-867).

Handling & Storage

Use/Stability	Stable for at least 6 months after receipt when stored at -20°C.
Handling	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -20°C.
Long Term Storage	-20°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
Concentration	1mg/ml (after reconstitution).
Formulation	Lyophilized. Contains PBS.
MW	104kDa (observed)
Purity	≥90% (SDS-PAGE)
Reconstitution	Reconstitute with 50µl sterile water. Further dilutions should be made with medium containing 5% fetal calf serum or a carrier protein.
Source	Produced in CHO cells. Human LRRC32 (aa 20-627) is fused to the Fc portion of human IgG1.
UniProt ID	Q14392

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)