# 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 ( $T_{req}$ ) cells.

## **Ordering Information**

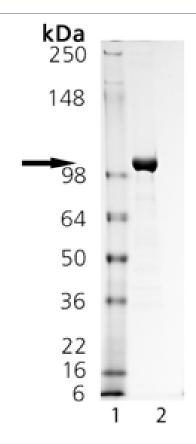
Order Online »

ALX-522-117-C050

50µg

Manuals, SDS & CofA

**View Online »** 



SDS-PAGE analysis: Lane 1: MW Marker, Lane 2: 2µg 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).

Endotoxin Content <0.1EU/µg purified protein (LAL test; Cape Cod

Associates).

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

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

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