

PGRP-S (human), (recombinant) (His- tag)

Peptidoglycan recognition protein 1, also known as Peptidoglycan recognition protein short, PGRP-S, PGLYRP1, PGLYRP, PGRP and TNFSF3L, is a secreted protein. It is highly expressed in bone marrow and weakly expressed in kidney, liver, small intestine, spleen, thymus, peripheral leukocyte, lung, fetal spleen and neutrophils. Peptidoglycan recognition proteins (PGRPs or PGLYRPs) are innate immunity proteins that are conserved from insects to mammals, recognize bacterial peptidoglycan, and function in antibacterial immunity and inflammation. Mammals have four PGRPs: PGLYRP1, PGLYRP2, PGLYRP3, and PGLYRP4. All PGRPs recognize bacterial peptidoglycan. The PGRPs likely play a role both in antibacterial defenses and several inflammatory diseases. They modulate local inflammatory responses in tissues (such as arthritic joints) and there is evidence for association of PGRPs with inflammatory diseases, such as psoriasis.

Ordering Information

[Order Online »](#)

ENZ-PRT322-0100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

kDa

M

116

66.2

45.0

35.0

25.0

18.4

14.4



As a result of glycosylation, the recombinant protein migrates as an approximately 40.5 kDa protein in SDS-PAGE under reducing conditions.

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name PGLYRP1, Peptidoglycan recognition protein short, TNFSF3L

Endotoxin Content <1 EU/μg (LAL method)

Formulation Lyophilized from sterile PBS, pH 7.4.

MW ~25.1 kDa (SDS-PAGE)

Purity ≥95% (SDS-PAGE)

Reconstitution Reconstitute with sterile deionized water. Reconstitution instructions are lot specific.

Source Produced in HEK293 cells. A DNA sequence encoding the human PGLYRP1 (Met1-Pro196) was expressed with a polyhistidine tag at the C-terminus.

UniProt ID O75594



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)