

GGT1 (human), (recombinant) (Fc-tag)

GGT1 belongs to the gamma-glutamyltransferase protein family. GGT1 catalyzes the transfer of the glutamyl moiety of glutathione to a variety of amino acids and dipeptide acceptors. GGT1 also initiates extracellular glutathione (GSH) breakdown, provides cells with a local cysteine supply and contributes to maintain intracellular GSH level. As part of the cell antioxidant defense mechanism, GGT1 can be detected in fetal and adult kidney and liver, adult pancreas, stomach, intestine, placenta and lung. Defects in GGT1 can cause glutathionuria which is known as an autosomal recessive disease.

Ordering Information

[Order Online »](#)

ENZ-PRT329-0100	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 132.5 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 CD224, Glutathione hydrolase 1

Endotoxin Content <1 EU/µg (LAL method)

Formulation Lyophilized from sterile PBS, pH 7.4.

MW ~122.2/59.5 kDa (SDS-PAGE)

Purity ≥90% (SDS-PAGE)

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

Source Produced in HEK293 cells. A DNA sequence encoding the human GGT1 (Pro27-Tyr569) was expressed with the fused Fc region of human IgG1 at the C-terminus.

UniProt ID P19440-1



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)