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 intracelular 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.

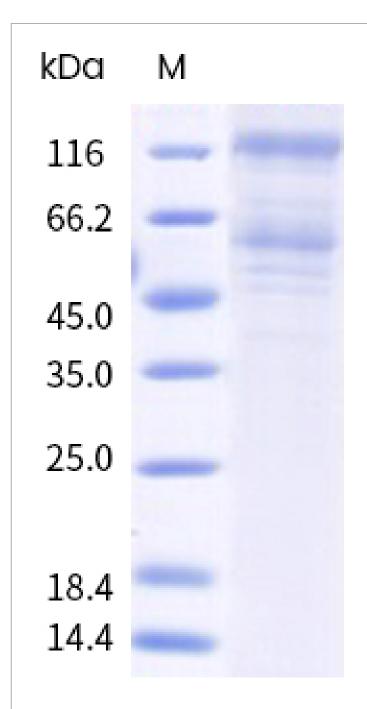
Ordering Information

Order Online »

ENZ-PRT329-0100 100µg

Manuals, SDS & CofA

View Online »



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

Last modified: November 7, 2024

