# GRO-β (human), (recombinant)

Growth regulated protein  $\beta$  (GRO- $\beta$ ), also known as CXCL2, is a chemokine that is secreted by macrophages, neutrophils, and monocytes. GRO- $\beta$  functions as a chemoattractant for leukocytes and hematopoietic stem cells. GRO- $\beta$  activity is mediated through binding the G-protein-coupled chemokine receptor CXCR2.

## **Ordering Information**

Order Online »

ENZ-PRT237-0010	10µg
ENZ-PRT237-0100	100µg

Manuals, SDS & CofA

**View Online** »

### **Handling & Storage**

**Handling** Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -80°C. For

long term storage, it is recommended to dilute to working aliquots in a 0.1% BSA

solution.

Long Term Storage -80°C

**Shipping** Ambient Temperature

## Regulatory Status RUO - Research Use Only

#### **Product Details**

Alternative Name MGSAβ, CXCL2, MIP-2α, GRO2

Endotoxin Content ≤1 EU/μg protein (LAL test)

Formulation Lyophilized from a sterile (0.2 micron) filtered solution

containing 0.1% Trifluoroacetic Acid (TFA).

**MW** ~7.9 kDa

Purity ≥95% (SDS-PAGE)

**Reconstitution** Reconstitute in sterile water at 0.1 mg/mL.

**Source** Produced in *E. coli*. Contains 73 amino acids.

UniProt ID P19875

Last modified: May 29, 2024



eu@enzolifesciences.com