

# GRO- $\beta$ (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	10 $\mu$ g
ENZ-PRT237-0100	100 $\mu$ g

## 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	MGSA $\beta$ , CXCL2, MIP-2 $\alpha$ , GRO2
Endotoxin Content	$\leq 1$ EU/ $\mu$ g protein (LAL test)
Formulation	Lyophilized from a sterile (0.2 micron) filtered solution containing 0.1% Trifluoroacetic Acid (TFA).
MW	~7.9 kDa
Purity	$\geq 95\%$ (SDS-PAGE)
Reconstitution	Reconstitute in sterile water at 0.1 mg/mL.
Source	Produced in <i>E. coli</i> . Contains 73 amino acids.
UniProt ID	P19875



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-  
fr@enzolifesciences.com](mailto:info-fr@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)