

# $\beta$ -Defensin 3 (human), (recombinant)

$\beta$ -Defensin 3 (BD-3) is a member of the defensin class of antimicrobial peptides that exert host defense responses against bacteria, fungi and viruses through microbial membranes. Interferon-gamma stimulates BD-3 expression and is important during adaptive immunity. BD-3 activates monocytes and mast cells, and has antibacterial functions towards Gram-negative and Gram-positive bacteria. Additionally, BD-3 blocks human immunodeficiency virus type 1 (HIV-1) replication through the downregulation of CXCR4, the HIV-1 co-receptor.

## Ordering Information

[Order Online »](#)

ENZ-PRT197-0020	20 $\mu$ g
-----------------	------------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	Reconstituted solution is stable for 3 months when stored at -20°C to -80°C and for 1 month when stored at 4°C after reconstituting as directed.
<b>Handling</b>	Centrifuge vial before opening. Suspend the product by gently pipetting the above recommended solution and allow several minutes for complete reconstitution. Do not vortex. For prolonged storage, dilute to working aliquots in a 0.1% BSA solution and store at -80°C. Avoid freeze/thaw cycles.
<b>Short Term Storage</b>	-20°C
<b>Long Term Storage</b>	-80°C
<b>Shipping</b>	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	BD-3
<b>Endotoxin Content</b>	≤1 EU/μg (LAL)
<b>Formulation</b>	Lyophilized from a sterile (0.2 micron) filtered aqueous solution containing 0.1% trifluoroacetic acid (TFA).
<b>MW</b>	5.2 kDa (with 45 amino acids)
<b>Purity</b>	≥95% (SDS-PAGE)
<b>Reconstitution</b>	Reconstitute with sterile 10 mM acetic acid at 0.1 mg/mL.
<b>Sequence</b>	GIINTLQKYYCRVRGGRC AVL SCLPKEEQIGKCSTRGRKCCRRKK
<b>Source</b>	Produced in <i>E. coli</i> .
<b>UniProt ID</b>	P81534



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)