# β-Defensin 3 (human), (recombinant)

β-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 Gramnegative 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

20µg

Manuals, SDS & CofA

**View Online »** 

## **Handling & Storage**

**Use/Stability** 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.

**Handling** 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.

Short Term Storage -20°C

Long Term Storage -80°C

**Shipping** Ambient Temperature

## Regulatory Status RUO - Research Use Only

### **Product Details**

Alternative Name BD-3

Endotoxin Content ≤1 EU/µg (LAL)

Formulation Lyophilized from a sterile (0.2 micron) filtered aqueous

solution containing 0.1% trifluoroacetic acid (TFA).

**MW** 5.2 kDa (with 45 amino acids)

Purity ≥95% (SDS-PAGE)

**Reconstitution** Reconstitute with sterile 10 mM acetic acid at 0.1 mg/mL.

Sequence GIINTLQKYYCRVRGGRCAVLSCLPKEEQIGKCSTRGRKCCRF

Source Produced in *E. coli*.

UniProt ID P81534



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com