## Defensins (human neutrophils)

A-defensins purified from human neutrophils. Defensins are short peptides secreted by neutrophils and epithelial cells as proproteins and activated by proteolysis. They have antibacterial, antiviral, and antifungal properties, and they promote immune response. Can be used in tissue culture or *in vivo*.

## **Ordering Information**

Order Online »

BML-SE514-0050

50µg

Manuals, SDS & CofA

View Online »

## **Handling & Storage**

**Use/Stability** As indicated on product label or CoA when stored as recommended.

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

## **Product Details**

Alternative Name HNPs, Human neutrophil peptides

MW 3-4

Purity ≥95%

Last modified: May 29, 2024

