



Product Data Sheet

Product Name: HNP-1, Defensin Human Neutrophil Peptide-1
Catalog Number: AS-60743 (0.5 mg) **Lot Number:** See label on vial
Sequence: H-Ala-Cys-Tyr-Cys-Arg-Ile-Pro-Ala-Cys-Ile-Ala-Gly-Glu-Arg-Arg-Tyr-Gly-Thr-Cys-Ile-Tyr-Gln-Gly-Arg-Leu-Trp-Ala-Phe-Cys-Cys-OH
(Disulfide bridge: 2-30, 4-19, 9-29) (3-letter code)
ACYCRIPACIAGERRYGTCTIYQGRLWAFCC
(Disulfide bridge: 2-30, 4-19, 9-29) (1-letter code)
Molecular Weight: 3443.1
% Peak Area by HPLC: ≥ 95
Appearance: Lyophilized white powder
Peptide Reconstitution: Using H₂O, reconstitute by adding 30-50 µl to 0.1 mg Defensin Human Neutrophil peptide.

Storage: Defensin Human Neutrophil peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at -20°C or lower. Reconstituted peptide can be aliquoted and stored at -20°C or lower.

Description: Mammalian defensins are abundant in the cytoplasmic azurophilic granules of neutrophils, Paneth cells of the small intestine and some macrophages. Human α-defensin-1 (HNP-1) is a peptide possessing both broad antimicrobial (both Gram-positive and Gram-negative bacteria) and cytotoxic activities. HNP-1 reduces adenoviral infection by more than 95%. Ref: Frick, I. et al. *J. Biol. Chem.* **278**, 16561 (2003); Valore, E. et al. *J. Clin. Invest.* **97**, 1624 (1996); Mizukawa, N. et al. *Anticancer Res.* **20**, 1125 (2000); Bastian, A. and H. Schafer, *Regul Pept.* **101**, 157 (2001).

Related Products:

Name	Cat #	Size
HNP-2, Defensin Human Neutrophil Peptide-2 (CYCRIPACIAGERRYGTCTIYQGRLWAFCC) (Disulfide bridge: 1-29, 3-18, 8-28)	AS-60744	0.1 mg
hBD-1, β-Defensin-1, human (DHYNCSVSSGGQCLYSACPIFTKIQGTCTYRGKAKCKK) (Disulfide bridge: 5-34, 12-27, 17-35)	AS-60740	0.1 mg
hBD-3, β-Defensin-3, human (GIINTLQKYCRVRRGRCVLSCLPKKEEQIGKCSTRGRKCCRKK) (Disulfide bridge: 11-40, 18-33, 23-41)	AS-60741	0.1 mg

For Research Use Only