



Product Data Sheet

Product Name:	hBD-1, β -Defensin-1, human	
Catalog Number:	AS-60740 (0.1 mg)	Lot Number: See label on vial
Sequence:	H-Asp-His-Tyr-Asn-Cys-Val-Ser-Ser-Gly-Gly-Gln-Cys-Leu-Tyr-Ser-Ala-Cys-Pro-Ile-Phe-Thr-Lys-Ile-Gln-Gly-Thr-Cys-Tyr-Arg-Gly-Lys-Ala-Lys-Cys-Cys-Lys-OH (Disulfide bridge: 5-34, 12-27, 17-35) DHYN CVSSGGQCLYSACPIFTKI QGTCYRGKAKCCK (Disulfide bridge: 5-34, 12-27, 17-35) (1-letter code)	
Molecular Weight:	3929.6	
Peptide Purity:	>95%	
Appearance:	Lyophilized white powder	

Peptide Reconstitution: hBD-1 peptide is freely soluble in water.

Storage: hBD-1 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.

Published Citations:

- Gryllos, I. et al. *PNAS* **105**, 16755 (2008).
- Kooi, C. & Sokol, PA. *Microbiol.* doi:10.1099/mic.0.028969-0 (2009).

Related Products:

Name	Cat #	Size
hBD-3, β -Defensin-3, human GIINTLQKYYCRVRGGRCVLSCLPKKEEQIGKCTRGRKCCRRKK (Disulfide bridge: 11-40, 18-33, 23-41)	AS-60741	0.1 mg
HNP-1, Defensin Human Neutrophil Peptide-1 ACYCRIPACIAGERRYGTCTIYQGRLWAFCC (Disulfide bridge: 2-30, 4-19, 9-29)	AS-60743	0.1 mg
HNP-2, Defensin Human Neutrophil Peptide-2 CYCRIPACIAGERRYGTCTIYQGRLWAFCC (Disulfide bridge: 1-29, 3-18, 8-28)	AS-60744	0.1 mg

For Research Use Only