

Product Data Sheet

| Product Name: | Elafin | | |
|--|---|-------------------------------|--|
| Catalog Number: | AS-61641 (0.1 mg) | Lot Number: See label on vial | |
| Sequence: | H-Ala-Gln-Glu-Pro-Val-Lys-Gly-Pro-Val-Ser-Thr-Lys-Pro-Gly-Ser-Cys- Pro-IIe-IIe-Leu-IIe-Arg-Cys-Ala-Met-Leu-Asn-Pro-Pro-Asn-Arg-Cys-Leu- Lys-Asp-Thr-Asp-Cys-Pro-Gly-IIe-Lys-Lys-Cys-Cys-Glu-Gly-Ser-Cys-Gly- Met-Ala-Cys-Phe-Val-Pro-Gln-OH (Disufide bonds between Cys16- Cys45, Cys23-Cys49, Cys32-Cys44, Cys38-Cys53) (3-letter code) AQEPVKGPVSTKPGSCPIILIRCAMLNPPNRCLKDTDCPGIKKCCEGSC GMACFVPQ (Disufide bonds between Cys16- Cys45, Cys23- Cys49, Cys32- Cys44, Cys38-Cys53) (1-letter code) | | |
| Molecular Weight: | 5999.3 | | |
| Peptide Purity: | >95% | | |
| Appearance: | Lyophilized white powder | | |
| Peptide Reconstitution: Elafin peptide is freely soluble in water. | | | |

Storage: PAR-2 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: This is a fragment of the elafin (elastase-specific inhibitor), an antiproteinase and antimicrobial (gram-positive and gram-negative respiratory pathogens) molecule that is expressed at epithelial sites. Elafin is a potent inhibitor of HNE and proteinase 3 produced in the skin, and in the airways, which is up-regulated in response to early inflammatory cytokines such as TNF and IL-1. Elafin, along with SLPI, also shares characteristics with antimicrobial defensinlike molecules in being a low molecular weight cationic peptide with the ability to eliminate pulmonary pathogens. Ref: King, A. et al. J. Clin. Endo. Metab. 88, 4426 (2003); Wiedow, O. et al . J. Biol. Chem. 265, 14791 (1990); Wiedow, O. et al. Biochem. Biophys. Res. Commun. 174, 6 (1991); Tremblay, G. et al. Am. J. Respir. Crit. Care Med. 154, 1092 (1996); Simpson, AJ. et al. J. Immunol. 167, 1778 (2001).

| Related Products: | | |
|----------------------------|----------|-------|
| Name | Cat # | Size |
| MeOSuc-Ala-Ala-Pro-Val-AFC | AS-24141 | 10 mg |
| MeOSuc-AAPV-AFC | | |

For Research Use Only