



Product Data Sheet

Product Name:	Endothelin 3, human, rat	
Catalog Number:	AS-24323 (0.5 mg) AS-24324 (1 mg)	Lot Number: See label on vial
Sequence:	H-Cys-Thr-Cys-Phe-Thr-Tyr-Lys-Asp-Lys-Glu-Cys-Val-Tyr-Tyr-Cys-His-Leu-Asp-Ile-Ile-Trp-OH (Disulfide bridge: 1-15 and 3-11) (3-letter code) CTCFITYKDKECVYYCHLDIIW (Disulfide bridge: 1-15 and 3-11) (1-letter code)	
Molecular Weight:	2643.1	
% Peak Area by HPLC:	≥ 95	
Appearance:	Lyophilized white powder	
Peptide Reconstitution:	Use distilled or higher quality water. Add water directly to the lyophilized peptide powder to obtain a final concentration of approximately 0.5 mg/mL to 1mg/mL or less. Gently vortex to mix. For peptides that have poor solubility in the suggested solvent, brief sonication may increase solubility in some cases.	
Storage:	Peptide is shipped at ambient temperature. Upon receipt, store lyophilized powder at –20°C or lower. Reconstituted peptide should be aliquoted into several freezer vials and stored at –20°C or lower. Do not freeze thaw.	

For Research Use Only