

Product Data Sheet

Product Name: Endothelin 3, human, rat

Catalog Number: AS-24323 (0.5 mg) Lot Number: See label on vial

AS-24324 (1 mg)

Sequence: H-Cys-Thr-Cys-Phe-Thr-Tyr-Lys-Asp-Lys-Glu-Cys-Val-Tyr-Tyr-Cys-His-

Leu-Asp-lle-lle-Trp-OH (Disulfide bridge: 1-15 and 3-11) (3-letter code) CTCFTYKDKECVYYCHLDIIW (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