



## Product Data Sheet

---

<b>Product Name:</b>	Endothelin 1, human, porcine	
<b>Catalog Number:</b>	AS-22859 (0.25 mg) AS-22860 (0.5 mg) AS-22861 (1 mg)	Lot Number: See label on vial
<b>Sequence:</b>	H-Cys-Ser-Cys-Ser-Ser-Leu-Met-Asp-Lys-Glu-Cys-Val-Tyr-Phe-Cys-His-Leu-Asp-Ile-Ile-Trp-OH (Disulfide bridge: 1-15 and 3-11) (3-letter code) CSCSSLMDKECVYFCHLDIIW (Disulfide bridge: 1-15 and 3-11) (1-letter code)	
<b>Molecular Weight:</b>	2492.0	
<b>Peptide Purity:</b>	>95%	
<b>Appearance:</b>	Lyophilized white powder	

**Peptide Reconstitution:** Initially try to dissolve a small amount of Endothelin peptide in water. If peptide does not dissolve, Endothelin peptide is freely soluble in 1% NH<sub>4</sub>OH.

**Storage:** Endothelin 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.

**Description:** ET-1 is a potent vasoconstrictor peptide derived from endothelial cells. It plays a role in regulation of cardiovascular functions. The binding of ET-1 to G-protein coupled receptor triggers a signal transduction cascade leading to activation of the protein kinase C (PKC). PKC inhibitor potentially reverses the up regulation of ET-1. The enhanced ET-1 expression associated with the activation of PKC occurs in early diabetes. Ref: Yanagisawa, M. et al. *Nature*. **332**, 411 (1988).

**Additional Information:** Listed below are relevant information that may provide a guideline on how to use this product. End users will have to adapt to their own specific applications.

Follicles (20–30 per vial) were treated with or without endothelin-1 (AnaSpec, San Jose, CA). To examine the role of cumulus cells in the endothelin-1-induced GVBD of oocytes, COCs were obtained by puncturing the largest follicles from preovulatory ovaries, whereas denuded oocytes were separated from cumulus cells by mechanical pipetting. These cells were cultured in minimum essential media (MEM, Invitrogen) without FBS in the presence of 4 mM hypoxanthine (Sigma) with or without 10 ng/ml endothelin-1 for 16 h at 37 °C in 5% CO<sub>2</sub>/95% air-[Kawamura, K. et al. \*Developmental Biol.\* \*\*327\*\*, 62 \(2009\).](#)

### Published Citations:

Ansar, S. et al. *J. Neurosurgery* **106**, DOI: 10.3171/jns.2007.106.5.881 (2006).  
Beg, SAS. et al. *J. Cer. Blood Flow Metab.* **26**, 846 (2006).

©AnaSpec, Inc. 34801 Campus Drive, Fremont, CA 94555  
Tel: (800)-452-5530 | [service@anaspec.com](mailto:service@anaspec.com) | [www.anaspec.com](http://www.anaspec.com)

Cervar-Zivkovic, M. et al. *Reprod. Sci.* **14**, 430 (2007).  
Wang, Y. et al. *Brain Res.* **1167**, 118 (2007).  
Kawamura, K. et al. *Developmental Biol.* **327**, 62 (2009).

**Related Products:**

<b>Name</b>	<b>Cat #</b>	<b>Size</b>
Endothelin 1, human, FAM-labeled FAM-CSCSSLMDKECVYFCHLDIIW (Disulfide bridge: 1-15 and 3-11)	AS-27116	0.5 mg
Big Endothelin-1 (1-38), human CSCSSLMDKECVYFCHLDIIWVNTPEHVVPYGLGSPRS (Disulfide bridge: 1-15 and 3-11)	AS-62946-05	0.5 mg
Big Endothelin 1 (1-39), porcine CSCSSLMDKECVYFCHLDIIWVNTPEHIVPYGLGSPRS (Disulfide bridge: 1-15 and 3-11)	AS-20715	1 mg

*For Research Use Only*