

# Big Endothelin-1 (human) (1-38)

Endothelin-converting enzyme substrate

Biological precursor of endothelin 1. Cleaved to yield endothelin-1 via the activity of the endothelin-converting enzyme (ECE). Endothelins are endothelium-derived vasoconstrictor peptides.

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-152-001-PC01	0.1mg*
------------------	--------

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Use/Stability** As indicated on product label or CoA when stored as recommended.

**Long Term Storage** -20°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Big ET-1 (1-38)

**Appearance** White to off-white powder.

**CAS** 121014-53-7

**Formula**  $C_{189}H_{282}N_{48}O_{56}S_5$

**MW** 4283.0

**Purity** ≥98% (HPLC)

**Quantity** \*Sold as net peptide weight.

**Sequence** H-Cys-Ser-Cys-Ser-Ser-Leu-Met-Asp-Lys-Glu-Cys-Val-Tyr-Phe-Cys-His-Leu-Asp-Ile-Ile-Trp-Val-Asn-Thr-Pro-Glu-His-Val-Val-Pro-Tyr-Gly-Leu-Gly-Ser-Pro-Arg-Ser-OH (Disulfide bonds between Cys<sup>1</sup>-Cys<sup>15</sup> and Cys<sup>3</sup>-Cys<sup>11</sup>)

**Solubility** 1mg/ml soluble in water; also soluble in 0.5% acetic acid.

**Last modified: May 29, 2024**



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)