

Endothelin-1 (human, mouse, rat, pig, bovine, dog)

Citations: 17

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-155-001-PC01	0.1mg*
ALX-155-001-PC05	0.5mg*
ALX-155-001-P001	1mg*

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Solution is stable for 1-2 weeks when stored at -20°C. Prepare solution just before use.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	ET-1
CAS	117399-94-7
Formula	$C_{109}H_{159}N_{25}O_{32}S_5$
Formulation	Lyophilized.
MI	14: 3576
MW	2492.0
Purity	≥98% (HPLC)
Quantity	*Sold as net peptide weight.
RTECS	KE9446500
Sequence	H-Cys-Ser-Cys-Ser-Ser-Leu-Met-Asp-Lys-Glu-Cys-Val-Tyr-Phe-Cys-His-Leu-Asp-Ile-Ile-Trp-OH (Disulfide bonds between Cys ¹ -Cys ¹⁵ and Cys ³ -Cys ¹¹)
Solubility	1mg/ml soluble in DMSO or 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-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

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