

# ET<sub>A</sub> receptor antagonist [BQ-610]

30-fold more potent ET<sub>A</sub> antagonist than BQ-123 in pig aortic smooth muscle membrane assay, same activity as BQ-123 when tested on pig cerebellum membrane.

## Ordering Information

[Order Online »](#)

|                  |      |
|------------------|------|
| ALX-155-011-P001 | 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

**CAS** 141595-53-1

**Formula**  $C_{36}H_{44}N_6O_6$

**MW** 656.8

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

**Sequence** (Hexahydro-1*H*-azepinyl)carbonyl-Leu-D-Trp(CHO)-D-Trp-OH

**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)