

# ET<sub>A</sub> receptor antagonist [BQ-123] . sodium salt

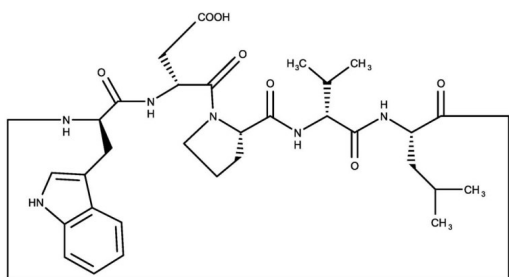
Highly potent and selective ET<sub>A</sub> antagonist.

Citations: 6

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Stable for at least 3 years after receipt when stored at -20°C. Stock solutions are stable for up to 3 months when stored at -20°C.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>CAS</b>	136553-81-6
<b>Formula</b>	$C_{31}H_{42}N_6O_7 \cdot Na$
<b>MW</b>	633.7
<b>Purity</b>	≥96%
<b>Quantity</b>	*Sold as net peptide weight.
<b>Sequence</b>	cyclo(D-Asp-Pro-D-Val-Leu-D-Trp . Na)
<b>Solubility</b>	Soluble in water.

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-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

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