

ET_A receptor antagonist [BQ-123] . sodium salt

Highly potent and selective ET_A antagonist.

Citations: 6

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-155-004-P001	1mg*
ALX-155-004-P005	5mg*

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	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.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

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

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-de@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-uk@enzolifesciences.com

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