

ET_B receptor antagonist . sodium salt

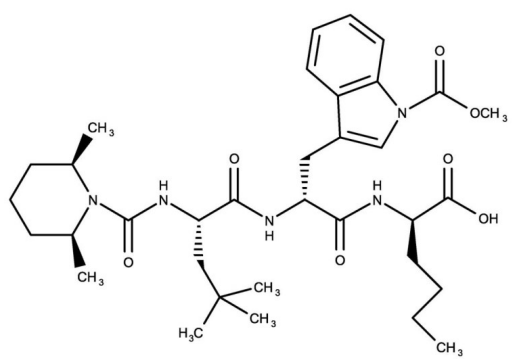
Highly potent and selective ET_B-receptor antagonist.

Citations: 2

[View Online »](#)

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	BQ-788 . Na
Appearance	White to off-white powder.
CAS	156161-89-6
Formula	$C_{34}H_{50}N_5O_7Na$
Formulation	Lyophilized from 0.1M sodium bicarbonate.
MW	663.7
Peptide Content	75-95%
Purity	≥95% (HPLC)
Sequence	N- <i>cis</i> -2,6-Dimethylpiperidinocarbonyl-L-γMeLeu-D-Trp(COOMe)-D-Nle-ONa
Solubility	Soluble in methanol (0.5mg/ml), DMSO (4mg/ml), 0.1M sodium bicarbonate (3mg/ml) or 100% ethanol (4mg/ml); insoluble 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-be@enzolifesciences.com

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

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

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