

Orexin A (human, bovine, rat, mouse)

Food intake modulator

Hypothalamic neuropeptide that stimulates food intake in rats with impressive potency. Binds specifically to the OX₁ receptor.

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-164-001-M001	1mg
ALX-164-0010	100µg

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

Formula $C_{152}H_{243}N_{47}O_{44}S_4$

Formulation Lyophilized.

MW 3561.2

Peptide Content 70-90%

Purity ≥98% (HPLC)

Sequence Glp-Pro-Leu-Pro-Asp-Cys-Cys-Arg-Gln-Lys-Thr-Cys-Ser-Cys-Arg-Leu-Tyr-Glu-Leu-Leu-His-Gly-Ala-Gly-Asn-His-Ala-Ala-Gly-Ile-Leu-Thr-Leu-NH₂
(Disulfide bonds between Cys⁶-Cys¹² and Cys⁷-Cys¹⁴).

Solubility Soluble in DMSO; can be diluted with buffers or distilled 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