

NBD peptide / NF-κB blocker (cell permeable)

For increased cell permeability the peptide was covalently linked to the 10 aa recognized by the TAT transporter.

Citations: 8

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-163-011-1	1Vial
---------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Keep on ice at all times. Protect from moisture. Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.
Short Term Storage	-20°C
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	NEMO-binding domain peptide
Concentration	10mM
Formula	$C_{121}H_{202}N_{48}O_{32}$
Formulation	Liquid. In sterile PBS.
MW	2841.2
Purity	≥95% (HPLC)
Quantity	50µl
Sequence	GRKKRRQRRR-GG-TALDWSWLQTE-amide

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)