

IKK-NBD peptide

NF-κB inhibitor

The peptide consists of two domains - an antennapedia homeodomain that confers cell permeability and the T735 to E745 region of IKKβ which is the NEMO binding domain (NBD). It is a selective inhibitor of NFκB activation. Blocks interaction of NEMO (an IκB complex regulatory protein) with the IκB kinase complex. The peptide (100 μM) inhibits E-selectin expression in human umbilical vein endothelial cells treated with TNFα and inhibits LPS-induced nitric acid release from a macrophage cell line. In an experimental mouse model of acute inflammation, IKK-NBD peptide (topically applied) inhibited PMA induced ear inflammation. Sold under license from Vanderbilt University

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-P607-0500	0.5mg
---------------	-------

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 Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Appearance Lyophilized solid

MW 3694.0

Purity ≥95% (HPLC)

Sequence Asp-Arg-Gln-Ile-Lys-Ile-Trp-Phe-Gln-Asn-Arg-Arg-Met-Lys-Trp-Lys-Lys-Thr-Ala-Leu-Asp-Trp-Ser-Trp-Leu-Gln-Thr-Glu

Solubility Soluble in DMSO (20mM). Note: This peptide will precipitate when added to tissue culture medium, even to 100µM. However, enough will remain in solution to enter many cell types.



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)