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 $\mathbf{C_{121}H_{202}N_{48}O_{32}}$

Formulation Liquid. In sterile PBS.

MW 2841.2

Purity ≥95% (HPLC)

Quantity 50μl

Sequence GRKKRRQRRR-GG-TALDWSWLQTE-amide

Last modified: May 29, 2024

