NF-κB SN50 cell permeable inhibitory peptide

Inhibits NF-kB translocation

Sequence contains the nuclear localization sequence (NLS residues 360-369) of the transcription factor NF- κ B p50 linked to a peptide cell-permeablization sequence, the hydrophobic region (h-region) of the signal peptide of Kaposi fibroblast growth factor (K-FGF). Peptide inhibits translocation of the NF- κ B active complex into the nucleus. In mouse endothelial LE-II cells induced with LPS, inhibition of NF- κ B nuclear translocation was 88% at 50 μ g/ml (18 μ M). In human monocytic THP-1 cells induced by LPS or TNF α there was 85% inhibition at 100 μ g/ml. Sold under license from Vanderbilt University.

Citations: 14

View Online »

Ordering Information

Order Online »

BML-P600-0005

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.

CAS 213546-53-3

MW 2781.5

Purity ≥97% (HPLC)

Sequence Ala-Ala-Val-Ala-Leu-Leu-Pro-Ala-Val-Leu-Leu-Ala-Leu-

Leu-Ala-Pro-Val-Gln-Arg-Lys-Arg-Gln-Lys-Leu-Met-Pro

Solubility Soluble in water, 100% ethanol or 50% ethanol.

Last modified: May 29, 2024



eu@enzolifesciences.com