

Product Data Sheet

Product Name: IKKy NEMO Binding Domain (NBD) Inhibitory Peptide

Catalog Number: AS-61169 (0.5 mg) Lot Number: See label on vial

Sequence: H-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-OH (3-letter code)

DRQIKIWFQNRRMKWKKTALDWSWLQTE (1-letter code)

Molecular Weight: 3693.3

Peptide Purity: >95%

Appearance: Lyophilized white powder

Peptide Reconstitution: IKKy NMB peptide is freely soluble in water.

Storage: IKKy NMB peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at -20°C or lower. Reconstituted peptide can be aliquoted and stored at -20°C or lower.

Description: A cell-permeable synthetic peptide NEMO-binding domain peptide (NBD peptide) corresponding to the NEMO amino-terminal alpha-helical region is shown to block TNF-alpha-induced NF-kB activation. The interaction of IKγNEMO with the IKK complex is critical for the activation of the IKK complex and the subsequent activation of NF-kB. Ref: Paria, B. et al. *J. Biol. Chem.* **278**, 37195 (2003); Tian, M. et al. *Blood* 10.1182/blood-2004-04-1248 (2005); Bharti, A. *Blood* **101**, 1053 (2003).

For Research Use Only