

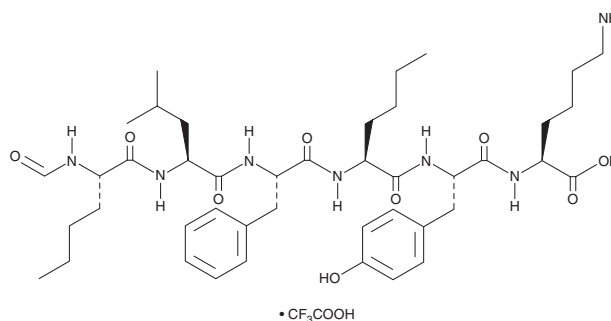
# PRODUCT INFORMATION



## N-formyl-Nle-Leu-Phe-Nle-Tyr-Lys (trifluoroacetate salt)

Item No. 35865

**CAS Registry No.:** 2918767-97-0  
**Formal Name:** N-formyl-L-norleucyl-L-leucyl-L-phenylalanyl-L-norleucyl-L-tyrosyl-L-lysine, monotrifluoroacetate salt  
**Synonym:** fNLFNYK  
**MF:** C<sub>43</sub>H<sub>65</sub>N<sub>7</sub>O<sub>9</sub> • CF<sub>3</sub>COOH  
**FW:** 938.1  
**Purity:** ≥98%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

N-formyl-Nle-Leu-Phe-Nle-Tyr-Lys (fNLFNYK) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the fNLFNYK (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. fNLFNYK (trifluoroacetate salt) is soluble in the organic solvent methanol.

### Description

fNLFNYK is a synthetic chemotactic peptide.<sup>1,2</sup> It binds to the formyl peptide receptor (FPR; K<sub>d</sub> = 2.2 nM).<sup>1</sup> fNLFNYK has been conjugated to various fluorophores to monitor FPR internalization.<sup>3</sup>

### References

1. Nidel, J.E. The formyl peptide chemotactic receptor: Cellular processing of peptide and receptor. *Receptor-Mediated Endocytosis. Receptors and Recognition*. Cuatrecasas, P. and Roth, T., editors, Springer (1983).
2. Boulay, F., Tardif, M., Brouchon, L., *et al.* Synthesis and use of a novel N-formyl peptide derivative to isolate a human N-formyl peptide receptor cDNA. *Biochem. Biophys. Res. Commun.* **168(3)**, 1103-1109 (1990).
3. Yan, P., Nanamori, M., Sun, M., *et al.* The immunosuppressant cyclosporin A antagonizes human formyl peptide receptor through inhibition of cognate ligand binding. *J. Immunol.* **177(10)**, 7050-7058 (2006).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 02/13/2024

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897

[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM