

# PRODUCT INFORMATION



## F1 (trifluoroacetate salt)

Item No. 35846

**Synonyms:** Tat-IKIP (46-60), Tat-Inhibitor of NF-κB  
Kinase-interacting Peptide

**Peptide Sequence:** YGRKKRRQRRRLSLAMTLGLAWLVF-OH      H-Tyr-Gly-Arg-Lys-Lys-Arg-Arg-Gln-Arg-Arg-Arg-Leu-Leu-Ser-Leu-Ala-Met-Thr-Leu-Gly-Leu-Ala-Trp-Leu-Val-Phe-OH

**MF:** C<sub>145</sub>H<sub>246</sub>N<sub>48</sub>O<sub>31</sub>S • XCF<sub>3</sub>COOH      • XCF<sub>3</sub>COOH

**FW:** 3,189.9

**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

F1 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the F1 (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

### Description

F1 is an anti-inflammatory peptide composed of the HIV-1 Tat protein transduction domain linked to a 15-amino acid peptide corresponding to residues 46-60 of inhibitor of NF-κB kinase-interacting peptide (IKIP).<sup>1</sup> It inhibits LPS-induced phosphorylation of IκB kinase α (IKKα) and IKKβ and nuclear translocation of NF-κB (p65) in isolated mouse peritoneal macrophages when used at a concentration of 5 μM. *In vivo*, F1 (5 mg/kg) decreases serum levels of IL-6, TNF-α, and IL-1β and increases survival in a mouse model of LPS-induced sepsis.

### Reference

1. Liu, H., Yan, Z., Zhao, Y., et al. A peptide derived from IKK-interacting protein attenuates NF-κB activation and inflammation. *J. Immunol.* 207(6), 1652-1661 (2021).

**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, 10/11/2023

### 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