

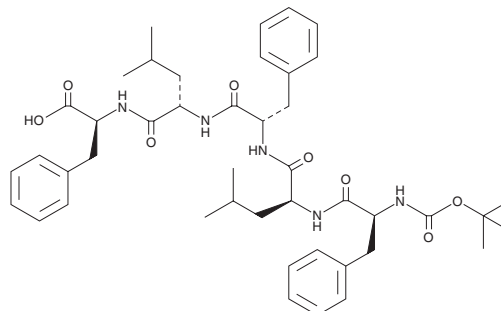
PRODUCT INFORMATION



Boc-FLFLF

Item No. 39673

CAS Registry No.: 73572-58-4
Formal Name: N-[(1,1-dimethylethoxy)carbonyl]-L-phenylalanyl-L-leucyl-L-phenylalanyl-L-leucyl-L-phenylalanine
Synonyms: Boc-Phe-Leu-Phe-Leu-Phe, Boc-PLPLP, L-BOC-2
Peptide Sequence: Boc-FLFLF-OH
MF: C₄₄H₅₉N₅O₈
FW: 786.0
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Boc-FLFLF is supplied as a solid. A stock solution may be made by dissolving the Boc-FLFLF in the solvent of choice, which should be purged with an inert gas. Boc-FLFLF is slightly soluble (0.1-1 mg/ml) in methanol and chloroform.

Description

Boc-FLFLF is a formyl peptide receptor (FPR) antagonist ($K_i = 1.46 \mu\text{M}$).¹ It inhibits increases in intracellular calcium levels, superoxide formation, and β -glucuronidase release induced by the FPR agonist N-formyl-Met-Leu-Phe (fMLP; Item No. 21495) in differentiated HL-60 cells (K_i s = 0.43, 1.04, and 1.82 μM , respectively).

Reference

1. Wenzel-Seifert, K., and Siefert, R. Cyclosporin H is a potent and selective formyl peptide receptor antagonist. Comparison with N-t-butoxycarbonyl-L-phenylalanyl-L-leucyl-L-phenylalanyl-L-leucyl-L-phenylalanine and cyclosporins A, B, C, D, and E. *J. Immunol.* **150**(10), 4591-4599 (1993).

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, 09/26/2024

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