

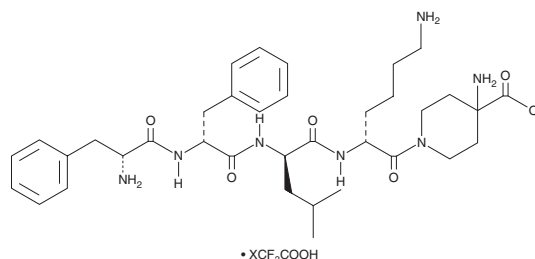
PRODUCT INFORMATION



Difelikefalin (trifluoroacetate salt)

Item No. 34542

CAS Registry No.: 2767426-81-1
Formal Name: N¹-(D-phenylalanyl-D-phenylalanyl-D-leucyl-D-lysyl)-4-amino-4-piperidinecarboxylic acid, trifluoroacetate salt
Synonyms: CR845, FE 202845
MF: C₃₆H₅₃N₇O₆ • XCF₃COOH
FW: 679.9
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

Difelikefalin (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the difelikefalin (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Difelikefalin (trifluoroacetate salt) is soluble in ethanol.

Description

Difelikefalin is a κ-opioid receptor (KOR) agonist.¹ It activates KOR in HEK293 cells expressing the human receptor (EC₅₀ = 0.16 nM in a transactivation assay) and inhibits forskolin-induced cAMP production in R1.G1 mouse thymoma cells (EC₅₀ = 0.048 nM). Difelikefalin is selective for KOR over the μ-opioid receptor (MOR; EC₅₀ = >1 μM in a transactivation assay). It reduces acetic acid-induced writhing, as well as scratching behavior induced by the KOR antagonist GNTI, in mice (ED₅₀s = 0.07 and 0.05 mg/kg, respectively).

Reference

1. Schteingart, C.D., Menzaghi, F., Jiang, G., *et al.* Synthetic peptide amides. *Cara Therapeutics, Inc.* US7,402,564B1 (2008).

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, 10/12/2023

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