

PRODUCT INFORMATION



Fz7-21 (trifluoroacetate salt)

Item No. 36922

Formal Name: N-acetyl-L-leucyl-L-prolyl-L-seryl-L- α -aspartyl-L- α -aspartyl-L-leucyl-L- α -glutamyl-L-phenylalanyl-L-tryptophyl-L-cysteinyl-L-histidyl-L-valyl-L-methionyl-L-tyrosynamide, trifluoroacetate salt

Synonym: dFz7-21

Peptide Sequence: Ac-LPSDDLEFWCHVMY-NH₂

MF: C₈₃H₁₁₄N₁₈O₂₃S₂ • XCF₃COOH

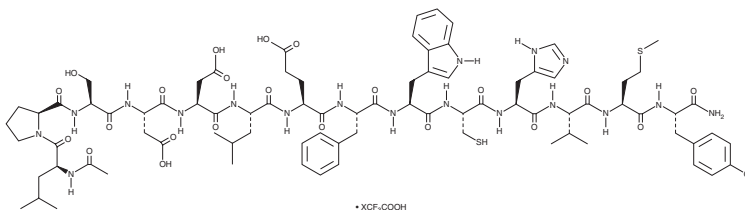
FW: 1796.0

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

Fz7-21 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the Fz7-21 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Fz7-21 (trifluoroacetate salt) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of Fz7-21 (trifluoroacetate salt) in these solvents is approximately 10 mg/ml.

Description

Fz7-21 is a peptide antagonist of the frizzled family G protein-coupled receptor 7 (FZD₇; IC₅₀ = 100 nM).¹ It inhibits Wnt3a-induced β -catenin signaling in a reporter assay using HEK293 cells. Fz7-21 (100 μ M) reduces the number of cells positive for leucine-rich α -2-glycoprotein 5 (Lrg5), a marker of intestinal stem cells, and the mRNA encoding achaete-scute homolog 2 (Ash-2) and Axin2 and increases mRNA encoding mucin 2 (Muc2) in primary mouse intestinal organoids.

Reference

1. Nile, A.H., de Sousa E. Melo, F., Mukund, S., *et al.* A selective peptide inhibitor of Frizzled 7 receptors disrupts intestinal stem cells. *Nat. Chem. Biol.* **14**(6), 582-590 (2018).

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, 11/16/2023

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