

PRODUCT INFORMATION



Fexapotide (trifluoroacetate)

Item No. 40716

CAS Registry No.: 1609252-56-3

Formal Name: L-isoleucyl-L- α -aspartyl-L-glutamyl-L-glutamyl-L-valyl-L-leucyl-L-seryl-L-arginyl-L-isoleucyl-L-lysyl-L-leucyl-L- α -glutamyl-L-isoleucyl-L-lysyl-L-arginyl-L-cysteinyl-L-leucine, triflutate

Synonym: NX-1207

Peptide Sequence: IDQQVLSRIKLEIKRCL-OH

MF: $C_{90}H_{163}N_{27}O_{25}S \cdot 5CF_3COOH$

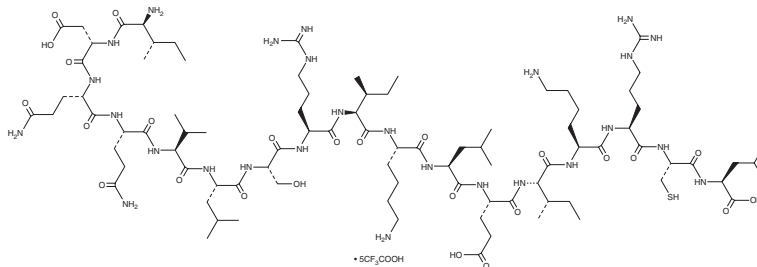
FW: 2,625.6

Purity: $\geq 95\%$

Supplied as: A solid

Storage: $-20^\circ C$

Stability: ≥ 4 years



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

Laboratory Procedures

Fexapotide (trifluoroacetate) is supplied as a solid. A stock solution may be made by dissolving the fexapotide (trifluoroacetate) in the solvent of choice, which should be purged with an inert gas. Fexapotide (trifluoroacetate) is soluble (≥ 10 mg/ml) in DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of fexapotide (trifluoroacetate) can be prepared by directly dissolving the solid in aqueous buffers. Fexapotide (trifluoroacetate) slightly soluble (0.1-1 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Fexapotide (trifluoroacetate) is a peptide.¹ It induces apoptosis in LNCaP prostate cancer cells when used at a concentration of 2.5 mg/ml. Intraprostatic administration of fexapotide triflutate (0.6 mg/animal per week) reduces the numbers of prostate glandular epithelial cells without affecting other cell types and decreases prostate volume in rats.

Reference

1. Averbach, P., Gohal, R., Fuska, M., *et al.* Fexapotide triflutate induces selective prostate glandular pharmaco-ablation in the rat. *Res. Rep. Urol.* **11**, 343-350 (2019).

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, 05/16/2024

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