

PRODUCT INFORMATION



FTSDVSKQMEEEEAVRLFIEWLKNGGPSSGAPPPS (trifluoroacetate salt)

Item No. 41332

Formal Name: (S)-N¹-((S)-6-amino-1-(((S)-3-hydroxy-1-(((3S,7S)-8-(((S)-3-hydroxy-1-(((2S,3R)-3-hydroxy-1-oxo-1-(((S)-1-oxo-3-phenylpropan-2-yl)amino)butan-2-yl)amino)-1-oxopropan-2-yl)amino)-7-(hydroxyamino)-2-methyl-4,5,8-trioxooctan-3-yl)amino)-1-oxopropan-2-yl)amino)-1-oxohexan-2-yl)-2-((S)-2-((4S,9S,14S,17S)-17-((S)-2-(2-((4S,7S)-7-((S)-2-((2S,3S)-2-((4S,7S)-7-((S)-2-((S)-6-amino-2-(2-((S)-1-amino-6-(2-((S)-1-(hydroxy-L-seryl-L-prolyl-L-prolyl-L-prolyl-L-alanylglycyl-L-seryl-L-seryl)pyrrolidine-2-carboxamido)acetamido)-1,4,5-trioxohexan-3-yl)hydrazineyl)hexanamido)-4-methylpentanamido)-4-(hydroxyamino)-8-(1H-indol-3-yl)-5,6-dioxooctanamido)-3-methylpentanamido)-3-phenylpropanamido)-1-((diaminomethylene)amino)-9-methyl-5,6-dioxodecan-4-yl)hydrazineyl)-3-methylbutanamido)-4,9,14-tris(hydroxyamino)-5,6,10,11,15,16-hexaoxooctadecanamido)-4-(methylthio)butanamido)pentanediamide, trifluoroacetate salt

MF: C₁₆₄H₂₅₂N₄₂O₅₃S • XCF₃COOH

FW: 3,692.1

Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

H—Phe—Thr—Ser—Asp—Val—Ser—Lys—Gln—Met—Glu—

Glu—Glu—Ala—Val—Arg—Leu—Phe—Ile—Glu—Trp—

Leu—Lys—Asn—Gly—Gly—Pro—Ser—Ser—Gly—Ala—

Pro—Pro—Pro—Ser—OH

• XCF₃COOH

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

Laboratory Procedures

FTSDVSKQMEEEEAVRLFIEWLKNGGPSSGAPPPS (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the FTSDVSKQMEEEEAVRLFIEWLKNGGPSSGAPPPS (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. FTSDVSKQMEEEEAVRLFIEWLKNGGPSSGAPPPS (trifluoroacetate salt) is soluble (≥10 mg/ml) in DMSO and sparingly soluble (1-10 mg/ml) in ethanol.

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 FTSDVSKQMEEEEAVRLFIEWLKNGGPSSGAPPPS (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. FTSDVSKQMEEEEAVRLFIEWLKNGGPSSGAPPPS (trifluoroacetate salt) is sparingly soluble (1-10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

FTSDVSKQMEEEEAVRLFIEWLKNGGPSSGAPPPS is a derivative of the endogenous incretin hormone glucagon-like peptide 1 (GLP-1; Item No. 24460).

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, 08/15/2024

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