

PRODUCT INFORMATION



Kisspeptin-234 (trifluoroacetate salt)

Item No. 35633

CAS Registry No.: 1848962-29-7

Formal Name: (S)-N¹-((6S,9R,15S)-9-((1H-indol-3-yl)methyl)-1-amino-6-(((S)-1-amino-1-oxo-3-phenylpropan-2-yl)carbamoyl)-15-benzyl-1-imino-8,11,14,17-tetraoxo-2,7,10,13,16-pentaazaocetadecan-18-yl)-2-((S)-2-((S)-2-((R)-2-acetamidopropanamido)-4-amino-4-oxobutanamido)-3-(1H-indol-3-yl)propanamido)succinamide, 2,2,2-trifluoroacetic acid

Ac—Ala—Asn—Trp—Asn—Gly—Phe—Gly—Trp—Arg—Phe—NH₂
• XCF₃COOH

Synonym: Kp-234

MF: C₆₃H₇₈N₁₈O₁₃ • XCF₃COOH

FW: 1,295.4

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

Kisspeptin-234 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the kisspeptin-234 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Kisspeptin-234 (trifluoroacetate salt) is slightly soluble in DMSO and dimethyl formamide.

Description

Kisspeptin-234 is a peptide antagonist of the kisspeptin receptor, previously known as G protein-coupled receptor 54 (GPR54).¹ It inhibits kisspeptin-10-induced inositol phosphate production in CHO cells expressing the human kisspeptin receptor (IC₅₀ = 7 nM) and kisspeptin-10-induced firing of isolated mouse gonadotropin-releasing hormone (GnRH) neurons. *In vivo*, kisspeptin-234 inhibits pulsatile GnRH secretion in pubertal female rhesus monkeys. It also inhibits kisspeptin-10-induced luteinizing hormone secretion in adult male rats and mice.

Reference

1. Roseweir, A.K., Kauffman, A.S., Smith, J.T., *et al.* Discovery of potent kisspeptin antagonists delineate physiological mechanisms of gonadotropin regulation. *J. Neurosci.* **29**(12), 3920-3929 (2009).

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, 04/18/2023

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