

PRODUCT INFORMATION



L-K6L9 (trifluoroacetate salt)

Item No. 38847

Formal Name: (S)-6-amino-N-((4S,7S,10S,13S,16S,19S,22S,25S,28S,31S)-35-amino-10,19,22-tris(4-aminobutyl)-4-carbamoyl-7,13,16,25,28-pentaisobutyl-2-methyl-6,9,12,15,18,21,24,27,30-nonaaxo-5,8,11,14,17,20,23,26,29-nonaazapentatriacontan-31-yl)-2-(((S)-2-(((S)-2-(((S)-6-amino-2-(((S)-2-amino-4-methylpentanamido)hexanamido)-4-methylpentanamido)-4-methylpentanamido)hexanamide, trifluoroacetate salt

Peptide Sequence: LKLLKLLKLLKLL-NH₂

MF: C₉₀H₁₇₄N₂₂O₁₅ • XCF₃COOH

FW: 1,804.5

Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

H—Leu—Lys—Leu—Leu—Lys—Lys—Leu—Leu—Lys—Lys—
Leu—Leu—Lys—Leu—Leu—NH₂
• XCF₃COOH

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

Laboratory Procedures

L-K6L9 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the L-K6L9 (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

L-K6L9 is a cytolytic peptide composed of L-isomers of leucine and lysine.¹ It is cytotoxic to LNCaP-CL1 androgen-independent, as well as 22Rv1 and LNCaP androgen-dependent, human prostate cancer cells (LC₅₀s = 4, 4, and 6 μM, respectively), as well as non-cancerous NIH3T3 mouse fibroblasts and OL human foreskin fibroblasts (LC₅₀s = 7 and 5 μM, respectively). It induces hemolysis of isolated human erythrocytes when used at a concentration of 100 μM.

Reference

1. Papo, N., Braunstein, A., Eshhar, Z., *et al.* Suppression of human prostate tumor growth in mice by a cytolytic D-, L-amino acid peptide: Membrane lysis, increased necrosis, and inhibition of prostate-specific antigen secretion. *Cancer Res.* **64**(16), 5779-5786 (2004).

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, 08/22/2023

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