

# PRODUCT INFORMATION



## ssK36 (trifluoroacetate salt)

Item No. 37514

**Formal Name:** (S)-2-(((S)-2-((6S,9S,18S, 21S,24S)-1- amino-6-(((S)- 2-((S)-3- amino-2- oxobutanoyl) pyrrolidin-1-yl)amino)-21- (4-amino butyl)-9- benzyl-24-(3-guanidino propyl)-1- imino-18- isopropyl-7, 10,13,16,19, 22,25-hepta oxo-2,8,11,14, 17,20,23-hepta azahexacosan- 26-oyl) pyrrolidin-1- yl) amino)-N<sup>1</sup>- ((6S,9S,12S)- 1,17-diamino- 9-(4-hydroxy benzyl)-6-(2- ((S)-1-hydroxy pyrrolidin- 2-yl)-2-oxo acetyl)-1,17- diimino-8,11- dioxo-2,7,10, 16-tetraaza heptadecan-12- yl)succinimide, trifluoroacetate salt

**Peptide Sequence:** APRFGGVKRPNRYRP-OH

**MF:** C<sub>79</sub>H<sub>127</sub>N<sub>29</sub>O<sub>18</sub> • XCF<sub>3</sub>COOH

**FW:** 1,771.1

**Purity:** ≥98%

**Supplied as:** A solid

**Storage:** -20°C

**Stability:** ≥4 years

H—Ala—Pro—Arg—Phe—Gly—Gly—Val—Lys—Arg—Pro—Asn—Arg—Tyr—Arg—Pro—OH  
• XCF<sub>3</sub>COOH

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

### Laboratory Procedures

ssK36 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the ssK36 (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

### Description

ssK36 is a peptide super-substrate for the histone methyltransferase SET domain-containing protein 2 (SETD2) corresponding to amino acids 29-43 of histone H3 lysine 36 (H3K36) with amino acid substitutions at positions 31, 32, 36, 37, and 39.<sup>1</sup> ssK36 is methylated 70-fold more and 290-fold faster than H3K36 (29-43) in cell-free assays.

### Reference

- Schuhmacher, M.K., Beldar, S., Khella, M.S., *et al.* Sequence specificity analysis of the SETD2 protein lysine methyltransferase and discovery of a SETD2 super-substrate. *Commun. Biol.* **3(1)**, 511 (2020).

#### 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, 10/11/2023

#### 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