

PRODUCT INFORMATION

Spike Glycoprotein (1147-1162) (SARS-CoV-2, SARS-CoV) (trifluoroacetate salt)

Item No. 37491

Formal Name: (3S,6S,9S,12S,15S,18S,21S,24S,27S,30,33S,36S)-9-((1H-imidazol-5-yl)methyl)-36-((S)-6-amino-2-((S)-2-((S)-2-amino-3-hydroxypropanamido)-3-phenylpropanamido)hexanamido)-12-(2-amino-2-oxoethyl)-15,24-bis(4-aminobutyl)-18-benzyl-33-(2-carboxyethyl)-27-(carboxymethyl)-21-(4-hydroxybenzyl)-6-((R)-1-hydroxyethyl)-3-(hydroxymethyl)-1-((S)-1-hydroxypyrrolidin-2-yl)-30-isobutyl-1,2,5,8,11,14,17,20,23,26,29,32,35-tridecaoxo-4,7,10,13,16,19,22,25,28,31,34-undecaazanonatriacontan-39-oic acid, trifluoroacetate salt

H—Ser—Phe—Lys—Glu—Glu—Leu—Asp—Lys—Tyr—Phe—
Lys—Asn—His—Thr—Ser—Pro—OH
• XCF₃COOH

Synonym: Surface Glycoprotein (1147-1162)

Peptide Sequence: SFKEELDKYFKNHTSP-OH

MF: C₉₀H₁₃₂N₂₂O₂₈ • XCF₃COOH

FW: 1,970.2

Purity: ≥95%

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

Spike glycoprotein (1147-1162) (SARS-CoV-2, SARS-CoV) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the spike glycoprotein (1147-1162) (SARS-CoV-2, SARS-CoV) (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

Spike glycoprotein (1147-1162) is a peptide corresponding to amino acids 1147-1162 of the severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) spike glycoprotein, also known as the surface glycoprotein.¹ This peptide is conserved across SARS-CoV, SARS-CoV-2, and various SARS-CoV-2 variants, including B.1.1.7, B.1.1.298, B.1.429, P.2, P.2, B.1.351, and B.1.617.2 and antibodies against this peptide neutralize SARS-CoV-2 and SARS-CoV pseudovirus particles in cellular assays.

Reference

- Li, T., Kan, Q., Ge, J., *et al.* A novel linear and broadly neutralizing peptide in the SARS-CoV-2 S2 protein for universal vaccine development. *Cell. Mol. Immunol.* **18**(11), 2563-2565 (2021).

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, 10/04/2023

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