

PRODUCT INFORMATION



SBP1 (trifluoroacetate salt)

Item No. 37469

Synonym: Spike Binding Peptide 1
Peptide Sequence: IEEQAKTFLDKFNHEAEDLFYQS-NH₂ H-Ile-Glu-Glu-Gln-Ala-Lys-Thr-Phe-Leu-Asp-
MF: C₁₂₇H₁₈₅N₃₁O₄₁ • XCF₃COOH Lys-Phe-Asn-His-Glu-Ala-Glu-Asp-Leu-Phe-
FW: 2,802.0 Tyr-Gln-Ser-NH₂
Purity: ≥95% • XCF₃COOH
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 binding peptide 1 (SBP1) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the SBP1 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. SBP1 (trifluoroacetate salt) is soluble in a 10% solution of acetonitrile/water.

Description

SBP1 is a peptide that corresponds to amino acids 21-43 of angiotensin-converting enzyme 2 (ACE2).¹ Lipid nanoparticles (LNPs) encapsulating the antiviral compound oseltamivir phosphate have been conjugated to SBP1 and studied *in vitro* to estimate their potential for providing controlled, long-term release of oseltamivir phosphate *in vivo*.¹ SBP1 (2%) immobilized to crosslinked networks of hydroxy acrylate and ethylxanthate ethyl acrylate, increases network capture efficiency of severe acute respiratory coronavirus 2 (SARS-CoV-2) spike glycoprotein, also known as surface glycoprotein, *in vitro*.²

References

1. Ucar, B., Acar, T., Arayici, P.P., *et al.* A nanotechnological approach in the current therapy of COVID-19: Model drug oseltamivir-phosphate loaded PLGA nanoparticles targeted with spike protein binder peptide of SARS-CoV-2. *Nanotechnology* **32(48)**, 485601 (2021).
2. Rahman, M.S., Rajawasam, C.W.H., De Alwis Watuthanthrige, N., *et al.* SARS-CoV-2 spike protein capture by peptide functionalized networks. *J. Polym. Sci.* **61(5)**, 391-397 (2023).

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, 10/13/2023

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