

PRODUCT INFORMATION



SP-A (196-215) (human) (trifluoroacetate salt)

Item No. 38838

Synonyms:	SPA4, Surfactant Protein A	
Peptide Sequence:	GDFRYSDGTPVNYTNWYRGE-OH	H—Gly—Asp—Phe—Arg—Tyr—Ser—Asp—Gly—Thr—Pro—
MF:	$C_{107}H_{145}N_{29}O_{35} \cdot XCF_3COOH$	
FW:	2,397.5	Val—Asn—Tyr—Thr—Asn—Trp—Tyr—Arg—Gly—Glu—OH
Purity:	≥98%	• XCF_3COOH
Supplied as:	A crystalline 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

SP-A (196-215) (human) (trifluoroacetate salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the SP-A (196-215) (human) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. SP-A (196-215) (human) (trifluoroacetate salt) is soluble in the organic solvent 10% acetonitrile/water.

Description

Surfactant protein A (SP-A) (196-215) is a synthetic peptide that corresponds to amino acids 196-215 of the C-terminal carbohydrate recognition domain of human SP-A.¹ It inhibits LPS-induced TNF- α release in JAWSII mouse dendritic cells when used at concentrations of 1 and 10 μ M. SP-A (196-215) (75 μ M) induces phagocytosis of *P. aeruginosa* by JAWSII cells.² Intratracheal administration of SP-A (196-215) (50 μ g/animal) decreases disease severity and lung colony forming units in a mouse model of *P. aeruginosa* infection.

References

1. Awasthi, S., Brown, K., King, C., *et al.* A toll-like receptor-4-interacting surfactant protein-A-derived peptide suppresses tumor necrosis factor- α release from mouse JAWS II dendritic cells. *J. Pharmacol. Exp. Ther.* **336**(3), 672-681 (2011).
2. Awasthi, S., Singh, B., Ramani, V., *et al.* TLR4-interacting SPA4 peptide improves host defense and alleviates tissue injury in a mouse model of *Pseudomonas aeruginosa* lung infection. *PLoS One* **14**(1), e0210979 (2019).

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, 08/01/2023

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