

PHC-SKKKK . 3 TFA

Negative control for Pam₃Cys-Ser-(Lys)₄ . trihydrochloride

Does not activate TLR2. Can be used as negative control for Pam3Cys-Ser-(Lys)₄ . trihydrochloride (Prod. No. ALX-165-066).

Lipopeptides are valuable tools for basic research in innate and acquired immunity. PHC-SKKKK (PHC-SK4) and Pam₃Cys-SKKKK (P3SK4) (Prod. No. ALX-165-066) have similar physicochemical properties, both have three fatty acid residues and the same peptide moiety. In contrast to P3SK4, PHCSK4 is not an activator of TLR2.

Citations: 4

[View Online »](#)

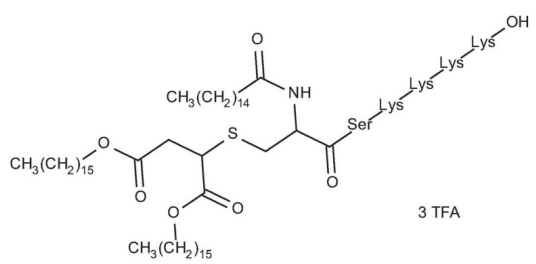
Ordering Information

[Order Online »](#)

ALX-165-071-M001	1mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -20°C.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	PHCSK4 . 3 TFA, N-palmitoyl-S-(1,2-bis-hexadecyloxy-carbonyl)-ethyl-(R)-cysteinyl-seryl-(lysyl)3-lysine . 3 CF3COOH
Appearance	Colorless powder
Concentration	1mg/ml after reconstitution.
Formula	C ₈₂ H ₁₅₈ N ₁₀ O ₁₃ S . 3 CF ₃ COOH
Formulation	Lyophilized.
MW	1524.3 . 342.1
Quality Control	Confirmed by ESI-MS.
Reconstitution	Reconstitute with 1 mL pyrogen-free deionized water in a pyrogen-free tube. Reconstitution can also be performed in DMSO (1mg/ml stock solution). Homogenize by vortexing or use a homogenizer or sonicator to prepare a homogenous solution. Further dilutions can be made with cell culture medium or pyrogen-free deionized water. PBS is not recommended for reconstitution or dilution.
Solubility	Soluble in pyrogen-free deionized water, DMSO or aqueous buffers at pH <7.5
Source	Synthetic



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)