

# PHC-SK4 . 3 TFA

Negative control for Pam<sub>3</sub>Cys-Ser-(Lys)<sub>4</sub> . trihydrochloride

Does not activate TLR2. Can be used as negative control for Pam<sub>3</sub>Cys-Ser-(Lys)<sub>4</sub> . trihydrochloride (Prod. No. ALX-165-066).

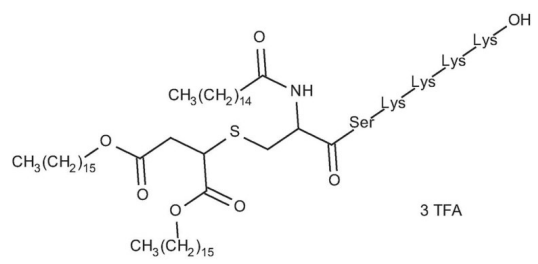
Lipopeptides are valuable tools for basic research in innate and acquired immunity. PHC-SK4 (PHC-SK4) and Pam<sub>3</sub>Cys-SK4 (P3SK4) (Prod. No. ALX-165-066) have similar physicochemical properties, both have three fatty acid residues and the same peptide moiety. In contrast to P3SK4, PHC-SK4 is not an activator of TLR2.

Citations: 4

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

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

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	PHCSK4 . 3 TFA, N-palmitoyl-S-(1,2-bis-hexadecyloxy-carbonyl)-ethyl-(R)-cysteinyl-seryl-(lysyl)3-lysine . 3 CF <sub>3</sub> COOH
<b>Appearance</b>	Colorless powder
<b>Concentration</b>	1mg/ml after reconstitution.
<b>Formula</b>	C <sub>82</sub> H <sub>158</sub> N <sub>10</sub> O <sub>13</sub> S . 3 CF <sub>3</sub> COOH
<b>Formulation</b>	Lyophilized.
<b>MW</b>	1524.3 . 342.1
<b>Quality Control</b>	Confirmed by ESI-MS.
<b>Reconstitution</b>	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.
<b>Solubility</b>	Soluble in pyrogen-free deionized water, DMSO or aqueous buffers at pH <7.5
<b>Source</b>	Synthetic

Last modified: May 29, 2024



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)