

Pam₃Cys-Ser-(Lys)₄ · trihydrochloride

Selective agonist of TLR1/TLR2 complex

Selective agonist of Toll-like receptor 1 (TLR1) complexed with TLR2. Cell permeable, water soluble synthetic cationic lipohexapeptide analog of the immunologically active N-terminal portion of bacterial lipoprotein that potentially activates monocytes and macrophages. Activates/represses an array of at least 140 genes partly involved in signal transduction and regulation of the immune response.

Citations: 31

[View Online »](#)

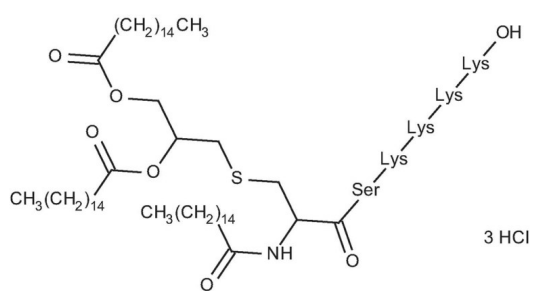
Ordering Information

[Order Online »](#)

ALX-165-066-M002	2mg
------------------	-----

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.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Pam3CSK4 . 3HCl, (S)-[2,3-Bis(palmitoyloxy)-(2-RS)-propyl]-N-palmitoyl-(R)-Cys-(S)-Ser-(S)-Lys4-OH . 3HCl, Pam3Cys-SKKKK
CAS	112208-00-1
Concentration	1mg/ml after reconstitution.
Formula	$C_{81}H_{156}N_{10}O_{13}S \cdot 3HCl$
Formulation	Lyophilized.
MW	1510.3 . 109.4
Quality Control	Confirmed by ESI-MS.
Reconstitution	Reconstitute with 2 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.
Technical Info / Product Notes	For negative control see Prod. No. ALX-165-071



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)