## 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 potently 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** 

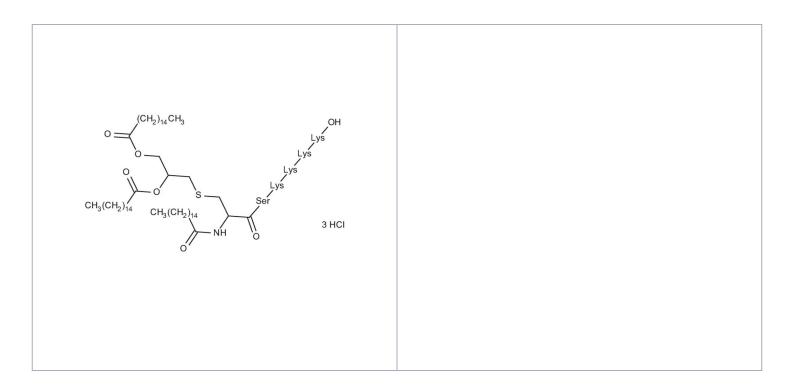
Order Online »

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  $C_{81}H_{156}N_{10}O_{13}S$  . 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

Last modified: May 29, 2024



info-

eu@enzolifesciences.com