

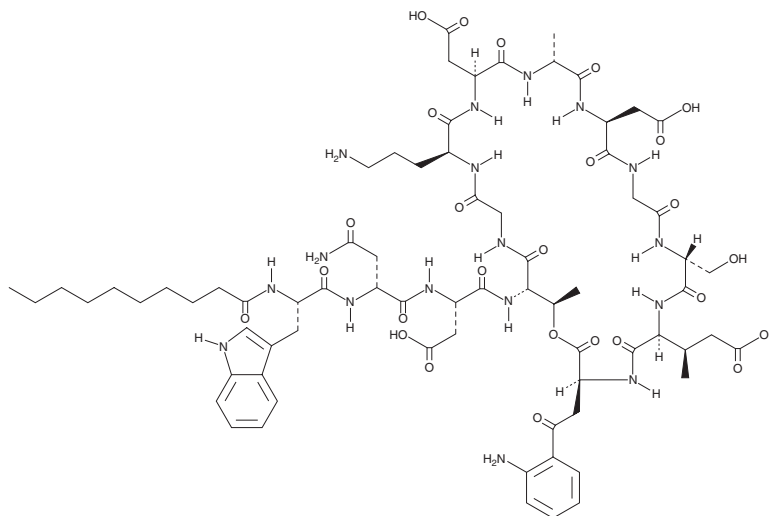
# PRODUCT INFORMATION



## Daptomycin Item No. 15615

**CAS Registry No.:** 103060-53-3  
**Formal Name:** 2,2'-((3S,6S,9R,15S,18R,21S,24S,30S,31R)-30-((S)-2-((R)-4-amino-2-((S)-2-decanamido-3-(1H-indol-3-yl)propanamido)-4-oxobutanamido)-3-carboxypropanamido)-3-(2-(2-aminophenyl)-2-oxoethyl)-24-(3-aminopropyl)-6-((R)-1-carboxypropan-2-yl)-9-(hydroxymethyl)-18,31-dimethyl-2,5,8,11,14,17,20,23,26,29-decaoxo-1-oxa-4,7,10,13,16,19,22,25,28-nonaazacyclohentriacontane-15,21-diyl) diacetic acid

**Synonym:** LY146032  
**MF:** C<sub>72</sub>H<sub>101</sub>N<sub>17</sub>O<sub>26</sub>  
**FW:** 1,620.7  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 223, 261, 370 nm  
**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

Daptomycin is supplied as a crystalline solid. A stock solution may be made by dissolving the daptomycin in the solvent of choice, which should be purged with an inert gas. Daptomycin is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of daptomycin in these solvents is approximately 5 and 1 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of daptomycin can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of daptomycin in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Daptomycin is a naturally occurring cyclic lipopeptide produced by the soil saprotroph *S. roseosporus* that is effective against resistant Gram-positive bacteria (MIC<sub>90</sub>s range from 0.25-16 mg/L).<sup>1,2</sup> It binds to bacterial plasma membranes to disrupt multiple aspects of function, including altering membrane potential and redirecting proteins essential for cell division and cell wall synthesis.<sup>2,3</sup>

### References

1. Tally, F.P. and DeBruin, M.F. *J. Antimicrob. Chemother.* **46**(4), 523-526 (2000).
2. Steenbergen, J.N., Alder, J., Throne, G.M., et al. *J. Antimicrob. Chemother.* **55**(3), 283-288 (2005).
3. Pogliano, J., Pogliano, N., and Silverman, J.A. *J. Bacteriol.* **194**(17), 4494-4504 (2012).

**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, 12/14/2022

### 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