

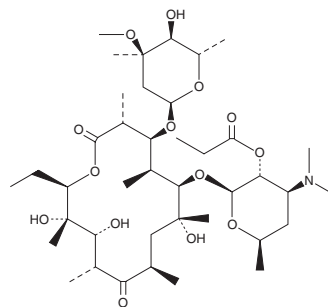
# PRODUCT INFORMATION



## Erythromycin 2'-Propionate

Item No. 31739

CAS Registry No.: 134-36-1  
Formal Name: 2'-propanoate erythromycin  
MF:  $C_{40}H_{71}NO_{14}$   
FW: 790.0  
Purity:  $\geq 95\%$   
Supplied as: A solid  
Storage:  $-20^{\circ}\text{C}$   
Stability:  $\geq 4$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Erythromycin 2'-propionate is supplied as a solid. A stock solution may be made by dissolving the erythromycin 2'-propionate in the solvent of choice, which should be purged with an inert gas. Erythromycin 2'-propionate is soluble in methanol and DMSO.

### Description

Erythromycin 2'-propionate is a macrolide antibiotic and an esterified form of erythromycin (Item No. 16486).<sup>1</sup> It is active against *S. aureus* when used at a concentration of 1  $\mu\text{g}/\text{ml}$ . Erythromycin 2'-propionate (1 mM) inhibits protein synthesis in a cell-free assay.

### Reference

1. Tardrew, P.L., Mao, J.C.H., and Kenney, D. Antibacterial activity of 2'-esters of erythromycin. *Appl. Microbiol.* **18(2)**, 159-165 (1969).

#### 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, 11/10/2022

#### 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