

# PRODUCT INFORMATION

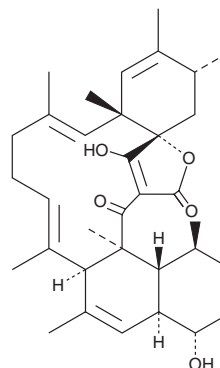


## Tetromycin B

Item No. 23766

**CAS Registry No.:** 180027-84-3  
**Formal Name:** (1S,4S,4aS,6aR,7E,11E,12aR,15R,16aS,20aS,20bR)-3,4,4a,6a,9,10,12a,15,16,20a,20b-dodecahydro-4,21-trihydroxy-1,6,7,11,12a,14,15,20a-octamethyl-8H-16a,19-metheno-16aH-benzo[b]naphth[2,1-j]oxacyclotetradecin-18,20(1H)-dione

**MF:** C<sub>36</sub>H<sub>46</sub>O<sub>5</sub>  
**FW:** 534.7  
**Purity:** ≥99%  
**Supplied as:** A 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

Tetromycin B is supplied as a solid. A stock solution may be made by dissolving the tetromycin B in the solvent of choice, which should be purged with an inert gas. Tetromycin B is soluble in organic solvents such as methanol, ethanol, DMSO, and dimethyl formamide.

### Description

Tetromycin B is a cysteine protease inhibitor with K<sub>i</sub> values of 0.62, 1.42, 32.5, and 1.59 μM for rhodesain, falcipain-2, cathepsin L, and cathepsin B, respectively.<sup>1</sup> It inhibits the growth of *T. brucei* *in vitro* (IC<sub>50</sub> = 30.87 μM). Tetromycin B is also cytotoxic to HEK293T kidney cells and J774.1 macrophages (IC<sub>50</sub>s = 71.77 and 20.2 μM, respectively).

### Reference

1. Pimentel-Elardo, S.M., Buback.V., Gulder, T.A.M., *et al.* New tetromycin derivatives with anti-trypanosomal and protease inhibitory activities. *Mar. Drugs* **9(10)**, 1682-1697 (2011).

**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, 09/30/2022

### 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