

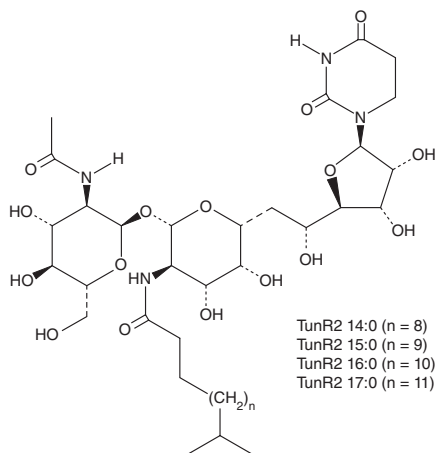
PRODUCT INFORMATION



TunR2

Item No. 31538

Formal Name: N-((2S,3R,4R,5R,6R)-2-(((2R,3R,4R,5S,6R)-3-acetamido-4,5-dihydroxy-6-(hydroxymethyl)tetrahydro-2H-pyran-2-yl)oxy)-6-((R)-2-((2R,3S,4R,5R)-5-(2,4-dioxotetrahydropyrimidin-1(2H)-yl)-3,4-dihydroxytetrahydrofuran-2-yl)-2-hydroxyethyl)-4,5-dihydroxytetrahydro-2H-pyran-3-yl)-12-methyltridecanamide
MF: $C_{39}H_{68}N_4O_{16}$ (for TunR2 16:0)
FW: 849.0
Purity: $\geq 95\%$ (mixture of congeners)
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.

Description

TunR2 is an antibiotic and derivative of tunicamycin (Item No. 11445).¹ It is active against *B. subtilis* (MIC = 0.3 $\mu\text{g/ml}$) and increases the efficacy of the β -lactam antibiotics oxacillin (Item No. 23954), methicillin (Item No. 21007), and penicillin G (Item No. 21615) against *B. subtilis* when used at a concentration of 0.4 $\mu\text{g/ml}$. Unlike tunicamycin, TunR2 is non-toxic to *S. cerevisiae* (MIC = $>10 \mu\text{g/ml}$) and does not inhibit glycosylation in a protein N-glycosylation assay. TunR2 also has reduced antiproliferative activity against MDA-MB-231 and CHO cells compared with tunicamycin.

Reference

1. Price, N.P., Hartman, T.M., Li, J., *et al.* Modified tunicamycins with reduced eukaryotic toxicity that enhance the antibacterial activity of β -lactams. *J. Antibiot. (Tokyo)* **70(11)**, 1070-1077 (2017).

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, 03/21/2024

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