

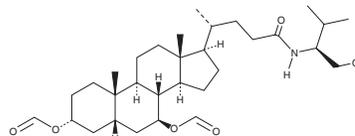
PRODUCT INFORMATION



A17

Item No. 39686

CAS Registry No.: 2568186-47-8
Formal Name: (3 α ,5 β ,7 β)-3,7-bis(formyloxy)-N-[(1S)-1-(hydroxymethyl)-2-methylpropyl]-cholan-24-amide
MF: C₃₁H₅₁NO₆
FW: 533.7
Purity: \geq 98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

A17 is supplied as a crystalline solid. A stock solution may be made by dissolving the A17 in the solvent of choice, which should be purged with an inert gas. A17 is sparingly soluble (1-10 mg/ml) in chloroform and slightly soluble (0.1-1 mg/ml) in methanol.

Description

A17 is a bile acid derivative.¹ It reduces triglyceride accumulation induced by oleic acid (Item Nos. 90260 | 24659), increases cell viability, and decreases oleic acid-induced lipid peroxidation in primary rat hepatocytes when used at concentrations of 10 and 25 μ M. A17 (25 μ M) reduces levels of CD36, also known as scavenger receptor B2 or fatty acid translocase, in primary rat hepatocytes. *In vivo*, A17 (50 mg/kg) reduces serum triglyceride, total cholesterol, and LDL levels, reverses hepatic steatosis, and decreases hepatic triglyceride, but not total cholesterol, levels in a hamster model of high-fat diet-induced non-alcoholic steatohepatitis (NASH).

Reference

1. Wang, Y., Zhu, Y., Niu, J., *et al.* A novel bile acid analog, A17, ameliorated non-alcoholic steatohepatitis in high-fat diet-fed hamsters. *Toxicol. Appl. Pharmacol.* **404**, 115169 (2020).

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, 04/04/2024

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