

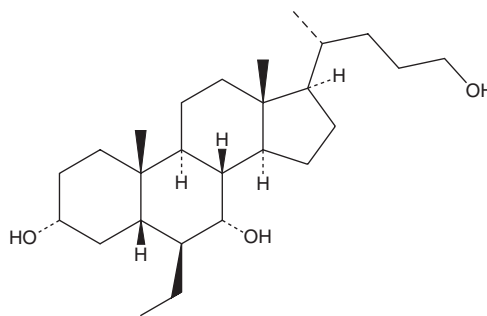
# PRODUCT INFORMATION



## BAR501 Impurity

Item No. 22461

**CAS Registry No.:** 1632118-70-7  
**Formal Name:** (3 $\alpha$ ,5 $\beta$ ,6 $\beta$ ,7 $\alpha$ )-6-ethyl-cholane-3,7,24-triol  
**MF:** C<sub>26</sub>H<sub>46</sub>O<sub>3</sub>  
**FW:** 406.6  
**Purity:**  $\geq$ 95%  
**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

BAR501 impurity is supplied as a crystalline solid. A stock solution may be made by dissolving the BAR501 impurity in the solvent of choice, which should be purged with an inert gas. BAR501 impurity is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of BAR501 impurity in ethanol and DMF is approximately 2 and 20 mg/ml, respectively. The solubility of BAR501 impurity in DMSO is approximately 100  $\mu$ g/ml.

BAR501 impurity is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BAR501 impurity should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. BAR501 impurity has a solubility of approximately 500  $\mu$ g/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

BAR501 impurity is an impurity found in the preparation of BAR501 (Item No. 22460) that acts as an agonist of the G protein-coupled bile acid-activated receptor (GP-BAR1). BAR501 impurity (10  $\mu$ M) induces a 150% increase in luciferase activity in a GP-BAR1 reporter gene assay.<sup>1</sup>

### Reference

1. Festa, C., Renga, B., D'Amore, C., *et al.* Exploitation of cholane scaffold for the discovery of potent and selective farnesoid X receptor (FXR) and G-protein coupled bile acid receptor 1 (GP-BAR1) ligands. *J. Med. Chem.* **57**(20), 8477-8495 (2014).

#### 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, 11/14/2022

#### 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