

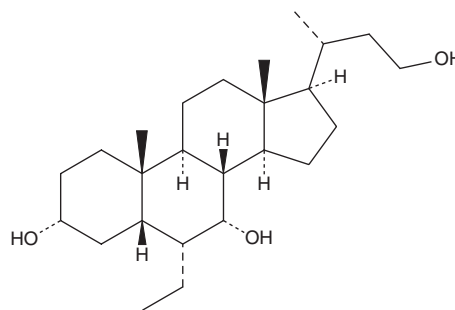
# PRODUCT INFORMATION



## BAR502

Item No. 22462

CAS Registry No.: 1612191-86-2  
Formal Name: (3 $\alpha$ ,5 $\beta$ ,6 $\alpha$ ,7 $\alpha$ )-6-ethyl-24-norcholeane-3,7,23-triol  
MF: C<sub>25</sub>H<sub>44</sub>O<sub>3</sub>  
FW: 392.6  
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

BAR502 is supplied as a crystalline solid. A stock solution may be made by dissolving the BAR502 in the solvent of choice, which should be purged with an inert gas. BAR502 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of BAR502 in these solvents is approximately 2, 0.1, and 20 mg/ml, respectively.

BAR502 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BAR502 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. BAR502 has a solubility of approximately 0.5 mg/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

BAR502 is a dual agonist of the G protein-coupled bile acid-activated receptor and farnesoid x receptor (GP-BAR1 and FXR; EC<sub>50</sub>s = 0.2 and 1  $\mu$ M, respectively).<sup>1</sup> It increases the expression of the FXR-regulated bile salt export pump (BSEP), organic solute transporter  $\alpha$  (OST $\alpha$ ), and small heterodimer partner (SHP) in primary hepatocytes isolated from FXR wild-type mice. BAR502 also increases cAMP-luciferase reporter gene expression, a marker of GP-BAR1 activity, in HEK293T cells.

### Reference

1. D'Amore, C., Di Leva, F.S., Sepe, V., *et al.* Design, synthesis, and biological evaluation of potent dual agonists of nuclear and membrane bile acid receptors. *J. Med. Chem.* **57**(3), 937-954 (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, 10/04/2022

#### 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