

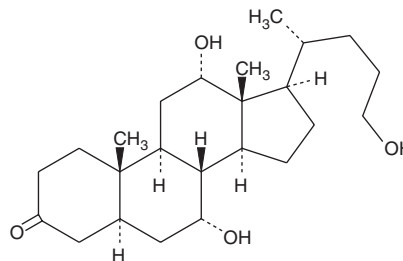
PRODUCT INFORMATION



3-keto Petromyzonol

Item No. 10007055

CAS Registry No.: 359436-56-9
Formal Name: 7 α ,12 α ,24-trihydroxy-5 α -cholane-3-one
MF: C₂₄H₄₀O₄
FW: 392.6
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

3-keto Petromyzonol is supplied as a crystalline solid. A stock solution may be made by dissolving the 3-keto petromyzonol in the solvent of choice, which should be purged with an inert gas. 3-keto Petromyzonol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 3-keto petromyzonol in these solvents is approximately 5, 14, and 20 mg/ml, respectively.

3-keto Petromyzonol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 3-keto petromyzonol should first be dissolved in DMF and then diluted with the aqueous buffer of choice. 3-keto Petromyzonol has a solubility of approximately 0.1 mg/ml in a 1:9 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

3-keto Petromyzonol is a synthetic intermediate useful for pharmaceutical synthesis.

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, 01/03/2023

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