

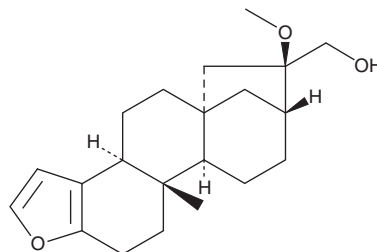
PRODUCT INFORMATION



16-O-Methylcafestol

Item No. 20259

CAS Registry No.: 108214-28-4
Formal Name: (3bS,5aS,7R,8R,10aR,10bS)-3b,4,5,6,7,8,9,10,10a,10b,11,12-dodecahydro-7-methoxy-10b-methyl-5a,8-methano-5aH-cyclohepta[5,6]naphtho[2,1-b]furan-7-methanol
MF: C₂₁H₃₀O₃
FW: 330.5
Purity: ≥98%
UV/Vis.: λ_{max}: 225 nm
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.

Laboratory Procedures

16-O-Methylcafestol is supplied as a crystalline solid. A stock solution may be made by dissolving the 16-O-methylcafestol in the solvent of choice. 16-O-Methylcafestol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of 16-O-methylcafestol in ethanol and DMSO is approximately 5 mg/ml and approximately 12 mg/ml in DMF.

16-O-Methylcafestol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 16-O-methylcafestol should first be dissolved in DMF and then diluted with the aqueous buffer of choice. 16-O-Methylcafestol has a solubility of approximately 0.02 mg/ml in a 1:50 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

16-O-Methylcafestol is a diterpene that has been found in robusta (*C. canephora*) coffee beans.¹ It has been used as a marker for robusta contamination and or adulteration of Arabica (*C. arabica*) roast coffee grounds in commercial production.

Reference

1. Gunning, Y., Defernez, M., Watson, A.D., *et al.* 16-O-methylcafestol is present in ground roast Arabica coffees: Implications for authenticity testing. *Food Chem.* **248**, 52-60 (2018).

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, 11/07/2022

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