

PRODUCT INFORMATION



MeTC7

Item No. 37338

CAS Registry No.: 1817841-22-7
Formal Name: 2-bromo-acetic acid, (6S,8aR,8bR,10aR,11R,13aR)-11-[(1R)-1,5-dimethylhexyl]-2,3,5,6,7,8,8a,8b,10,10a,11,12,13,13a-tetradecahydro-2,8a,10a-trimethyl-1,3-dioxo-4a,13b-etheno-1H,9H-benzo[c]cyclopenta[h][1,2,4]triazolo[1,2-a]cinnolin-6-yl ester

MF: C₃₂H₄₈BrN₃O₄

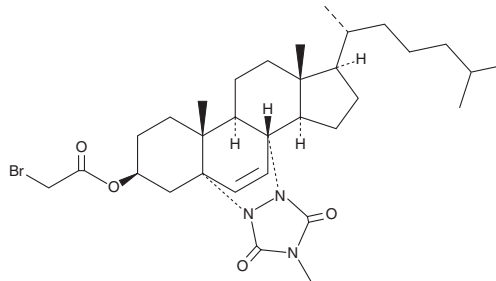
FW: 618.7

Purity: ≥95%

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥2 years



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

Laboratory Procedures

MeTC7 is supplied as a crystalline solid. A stock solution may be made by dissolving the MeTC7 in the solvent of choice, which should be purged with an inert gas. MeTC7 is soluble in the organic solvent ethanol at a concentration of approximately 1 mg/ml.

Description

MeTC7 is a vitamin D receptor antagonist ($IC_{50} = 2.9 \mu M$ in a fluorescence polarization assay).¹ It inhibits vitamin D receptor transactivation in HEK293 cells ($IC_{50} = 20.8 \mu M$). It decreases the viability of OVCAR-8, OV-2008, SKOV3, Caov-3, and IGROV-1 ovarian cancer cells in a concentration-dependent manner. MeTC7 (250 nM) reduces retinoid X receptor α (RXR α) and importin-4 protein levels in OV-2008 cells. It reduces tumor growth in an SH-SY5Y neuroblastoma mouse xenograft model when administered at a dose of 10 mg/kg.

Reference

1. Khazan, N., Kim, K.K., Hansen, J.N., *et al.* Identification of a vitamin-D receptor antagonist, MeTC7, which inhibits the growth of xenograft and transgenic tumors *in vivo*. *J. Med. Chem.* **65**(8), 6039-6055 (2022).

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, 09/29/2022

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