

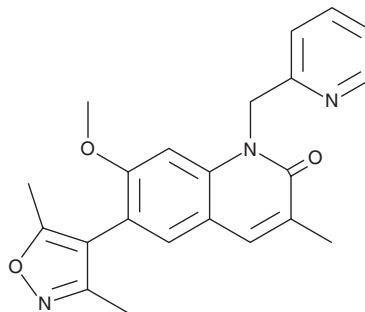
PRODUCT INFORMATION



ODM-207

Item No. 37757

CAS Registry No.: 1801503-93-4
Formal Name: 6-(3,5-dimethyl-4-isoxazolyl)-7-methoxy-3-methyl-1-(2-pyridinylmethyl)-2(1H)-quinolinone
MF: C₂₂H₂₁N₃O₃
FW: 375.4
Purity: ≥98%
UV/Vis.: λ_{max}: 213, 235, 331, 344 nm
Supplied as: A 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

ODM-207 is supplied as a solid. A stock solution may be made by dissolving the ODM-207 in the solvent of choice, which should be purged with an inert gas. ODM-207 is soluble in the organic solvent DMSO at a concentration of approximately 10 mM.

Description

ODM-207 is an inhibitor of bromodomain and extra terminal domain (BET) bromodomains. It reduces the proliferation of isolated human metastatic apocrine sweat gland cancer cells (IC₅₀ = 1.99 μM).¹ ODM-207, in combination with the poly(ADP-ribose) polymerase (PARP) inhibitor olaparib (Item No. 10621), decreases proliferation of, and induces DNA double-strand breaks in, isolated human metastatic apocrine sweat gland cancer cells deficient in partner and localizer of BRCA2 (PALB2; IC₅₀ = 650 nM).

Reference

1. Mäkelä, R., Härmä, V., Fajardo, N.B., *et al.* Ex vivo analysis of DNA repair targeting in extreme rare cutaneous apocrine sweat gland carcinoma. *Oncotarget* **12(11)**, 1100-1109 (2021).

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, 05/01/2023

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