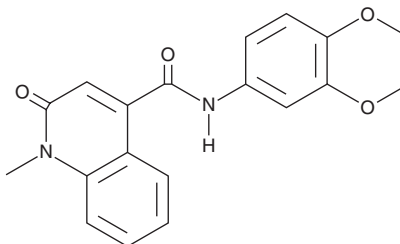


PRODUCT INFORMATION



CeMMEC13 Item No. 20224

CAS Registry No.: 1790895-25-8
Formal Name: N-(2,3-dihydro-1,4-benzodioxin-6-yl)-1,2-dihydro-1-methyl-2-oxo-4-quinolinecarboxamide
MF: C₁₉H₁₆N₂O₄
FW: 336.3
Purity: ≥98%
UV/Vis.: λ_{max}: 211, 337 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

CeMMEC13 is supplied as a crystalline solid. A stock solution may be made by dissolving the CeMMEC13 in the solvent of choice, which should be purged with an inert gas. CeMMEC13 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of CeMMEC13 in these solvents is approximately 20 mg/ml.

CeMMEC13 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CeMMEC13 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CeMMEC13 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CeMMEC13 is an isoquinolinone that selectively inhibits the second bromodomain of TAF1 (IC₅₀ = 2.1 μM).¹ It does not bind to bromodomains of BRD4, BRD9, or CREBBP.¹ CeMMEC13 synergizes with (+)-JQ1 (Item No. 11187) to inhibit the proliferation of THP-1 and H23 lung adenocarcinoma cells.¹

Reference

1. Sdelci, S., Lardeau, C.-H., Tallant, C., *et al.* Mapping the chemical chromatin reactivation landscape identifies BRD4-TAF1 cross-talk. *Nat. Chem. Biol.* **12**(7), 504-510 (2016).

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, 03/27/2025

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