

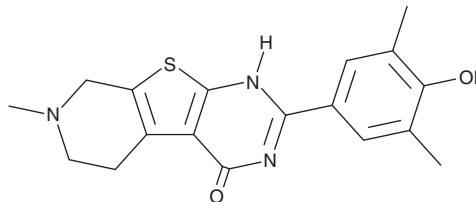
PRODUCT INFORMATION



FL-411

Item No. 37547

CAS Registry No.: 2118944-88-8
Formal Name: 5,6,7,8-tetrahydro-2-(4-hydroxy-3,5-dimethylphenyl)-7-methyl-pyrido[4',3':4,5]thieno[2,3-d]pyrimidin-4(1H)-one
Synonyms: BRD4-IN-1, Bromodomain-containing Protein 4 Inhibitor 1
MF: C₁₈H₁₉N₃O₂S
FW: 341.4
Purity: ≥98%
UV/Vis.: λ_{max}: 341.4
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

FL-411 is supplied as a solid. A stock solution may be made by dissolving the FL-411 in the solvent of choice, which should be purged with an inert gas. FL-411 is soluble in DMSO.

Description

FL-411 is an inhibitor of bromodomain-containing protein 4 (BRD4; IC₅₀ = 0.43 μM).¹ It is selective for BRD4 over BRD2 and BRD3 (IC₅₀s = 24.6 and >100 μM, respectively). FL-411 inhibits the proliferation of MCF-7 and MDA-MB-231 breast cancer cells (IC₅₀s = 5.48 and 8.29 μM, respectively) and induces autophagy in the same cells when used at concentrations of 3, 10, and 150 μM.

Reference

1. Ouyang, L., Zhang, L., Liu, J., *et al.* Discovery of a small-molecule bromodomain-containing protein 4 (BRD4) inhibitor that induces AMP-activated protein kinase-modulated autophagy-associated cell death in breast cancer. *J. Med. Chem.* **60(24)**, 9990-10012 (2017).

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, 04/06/2023

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