

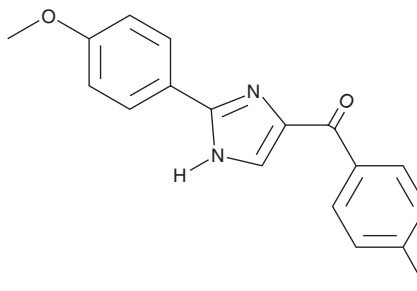
PRODUCT INFORMATION



CAY10647

Item No. 10791

CAS Registry No.: 1253697-49-2
Formal Name: (4-fluorophenyl)(2-(4-methoxyphenyl)-1H-imidazol-4-yl)methanone
MF: C₁₇H₁₃FN₂O₂
FW: 296.3
Purity: ≥98%
UV/Vis.: λ_{max}: 258, 325 nm
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

CAY10647 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10647 in the solvent of choice, which should be purged with an inert gas. CAY10647 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of CAY10647 in these solvents is approximately 3 mg/ml.

Description

CAY10647 is an aryl-benzoyl-imidazole which prevents proliferation of multidrug resistant cancer cells *in vitro* (IC₅₀ = 28-63 nM) and inhibits melanoma xenograph tumor growth *in vivo* (10-30 mg/kg).¹ This compound also binds to tubulin (K_i = 3.4 μM) and prevents tubulin polymerization.¹

Reference

1. Chen, J., Wang, Z., Li, C.M., *et al.* Discovery of novel 2-aryl-4-benzoyl-imidazoles targeting the colchicines binding site in tubulin as potential anticancer agents. *J. Med. Chem.* **53**(20), 7414-27 (2010).

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, 05/15/2023

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