

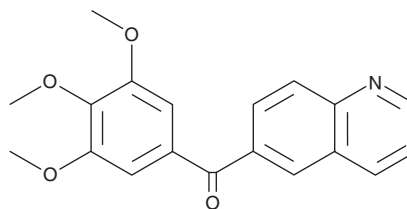
# PRODUCT INFORMATION



## MPT0B014

Item No. 21718

**CAS Registry No.:** 1215208-59-5  
**Formal Name:** 6-quinolinyl(3,4,5-trimethoxyphenyl)methanone  
**MF:** C<sub>19</sub>H<sub>17</sub>NO<sub>4</sub>  
**FW:** 323.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 250 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

MPT0B014 is supplied as a crystalline solid. A stock solution may be made by dissolving the MPT0B014 in the solvent of choice, which should be purged with an inert gas. MPT0B014 is soluble in the organic solvent DMSO at a concentration of approximately 5 mg/ml.

### Description

MPT0B014 is an inhibitor of tubulin polymerization.<sup>1</sup> It inhibits proliferation of A549, H1299, and H226 human non-small cell lung cancer (NSCLC) cells (IC<sub>50</sub>s = 0.041, 0.046, and 0.041 μM, respectively), as well as adriamycin-resistant NCI/ADR cells that overexpress P-glycoprotein (IC<sub>50</sub> = ~0.03 μM). MPT0B014 (0.3 μM) induces cell cycle arrest at the G<sub>2</sub>/M phase and apoptosis in A549 cells. It reduces tumor growth in an A549 mouse xenograft model when administered in combination with the EGFR tyrosine kinase inhibitor erlotinib (Item No. 10483).

### Reference

1. Tsai, A.-C., Pai, H.-C., Wang, C.-Y., *et al.* *In vitro* and *in vivo* anti-tumour effects of MPT0B014, a novel derivative aroylquinoline, and in combination with erlotinib in human non-small-cell lung cancer cells. *Br. J. Pharmacol.* **171**(1), 122-133 (2014).

#### 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, 12/16/2022

#### 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