

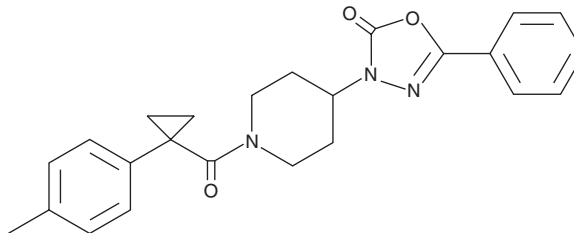
# PRODUCT INFORMATION



## ML-191

Item No. 15182

**CAS Registry No.:** 931695-79-3  
**Formal Name:** 5-phenyl-3-(1-(1-(*p*-tolyl)cyclopropane-1-carbonyl)piperidin-4-yl)-1,3,4-oxadiazol-2(3H)-one  
**Synonyms:** CCG-152883, CID-23612552  
**MF:** C<sub>24</sub>H<sub>25</sub>N<sub>3</sub>O<sub>3</sub>  
**FW:** 403.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 269 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

ML-191 is supplied as a crystalline solid. A stock solution may be made by dissolving the ML-191 in the solvent of choice, which should be purged with an inert gas. ML-191 is soluble in the organic solvent DMSO at a concentration of approximately 5 mg/ml.

### Description

ML-191 is an antagonist of GPR55.<sup>1</sup> It inhibits GPR55 signaling induced by lysophosphatidylinositol (LPI; EC<sub>50</sub> = 1.076 μM in U2OS cells overexpressing GPR55). ML-191 inhibits LPI-induced phosphorylation of ERK1/2 (IC<sub>50</sub> = 328 nM) and receptor-dependent translocation of PKCβII when used at a concentration of 30 μM.<sup>1</sup>

### Reference

1. Heynen-Genel, S., Dahl, R., Shi, S., *et al.* Selective GPR55 antagonists: Screening for selective ligands for GPR55. *Probe Reports from the NIH Molecular Libraries Program* (2011).

#### 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, 10/21/2022

#### 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