

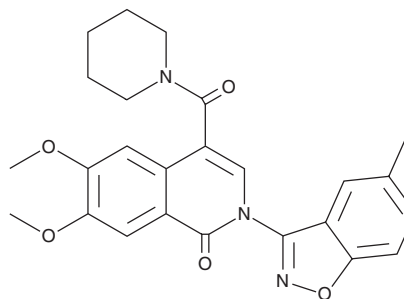
# PRODUCT INFORMATION



**AS-2717638**

Item No. 37787

**CAS Registry No.:** 2148339-28-8  
**Formal Name:** 6,7-dimethoxy-2-(5-methyl-1,2-benzisoxazol-3-yl)-4-(1-piperidinylcarbonyl)-1(2H)-isoquinolinone  
**MF:** C<sub>25</sub>H<sub>25</sub>N<sub>3</sub>O<sub>5</sub>  
**FW:** 447.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 253, 295 nm  
**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

AS-2717638 is supplied as a solid. A stock solution may be made by dissolving the AS-2717638 in the solvent of choice, which should be purged with an inert gas. AS-2717638 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of AS-2717638 in these solvents is approximately 2 mg/ml.

## Description

AS-2717638 is a lysophosphatidic acid receptor 5 (LPA<sub>5</sub>) antagonist (IC<sub>50</sub> = 0.038 μM).<sup>1</sup> It is selective for LPA<sub>5</sub> over LPA<sub>1</sub>, LPA<sub>2</sub>, and LPA<sub>3</sub> (IC<sub>50</sub>s = >10 μM for all) but also binds to adenosine A<sub>1</sub> and μ-opioid receptors at 10 μM. It induces analgesia in a rat model of allodynia induced by LPA or geranylgeranyl pyrophosphate (GGPP; Item No. 63330) when administered at doses of 3, 10, and 30 mg/kg. AS-2717638 (10 and 30 mg/kg) inhibits allodynia induced by prostaglandin E<sub>2</sub> (PGE<sub>2</sub>; Item No. 14010), PGF<sub>2α</sub>, or AMPA in rats. It increases the paw withdrawal threshold and latency in the von Frey hair and plantar tests, respectively, in rats when administered at a dose of 10 mg/kg.

## Reference

1. Murai, N., Hiyama, H., Kiso, T., *et al.* Analgesic effects of novel lysophosphatidic acid receptor 5 antagonist AS2717638 in rodents. *Neuropharmacology* **126**, 97-107 (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, 10/24/2022

## 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