

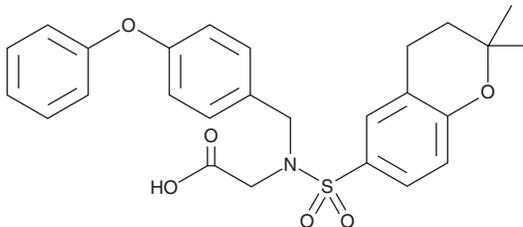
# PRODUCT INFORMATION



## LEI-106

Item No. 16419

**CAS Registry No.:** 1620582-23-1  
**Formal Name:** N-[(3,4-dihydro-2,2-dimethyl-2H-1-benzopyran-6-yl)sulfonyl]-N-[(4-phenoxyphenyl)methyl]-glycine  
**MF:** C<sub>26</sub>H<sub>27</sub>NO<sub>6</sub>S  
**FW:** 481.6  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 245 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

LEI-106 is supplied as a crystalline solid. A stock solution may be made by dissolving the LEI-106 in the solvent of choice, which should be purged with an inert gas. LEI-106 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of LEI-106 in these solvents is approximately 0.1, 30, and 50 mg, respectively.

LEI-106 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, LEI-106 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. LEI-106 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

In humans, two forms of diacylglycerol lipase, DAGL $\alpha$  and DAGL $\beta$ , generate the endocannabinoid 2-arachidonoyl glycerol (2-AG; Item No. 62160) by attacking DAG at the *sn*-1 position. LEI-106 is a potent *in vitro* inhibitor of *sn*-1 DAGL $\alpha$  (IC<sub>50</sub> = 18 nM).<sup>1</sup> It blocks the hydrolysis of *sn*-1-oleoyl-2-AG, the natural substrate of DAGL $\alpha$ , with a K<sub>i</sub> value of 0.7  $\mu$ M.<sup>1</sup> LEI-106 inhibits the hydrolysis of 2-AG by the monoacylglycerol lipase ABHD6 in mouse brain membrane homogenates and in HEK293T cell membrane preparations (K<sub>i</sub> = 0.8  $\mu$ M).<sup>1</sup>

### Reference

1. Janssen, F.J., Deng, H., Allarà, M., *et al.* Discovery of glycine sulfonamides as dual inhibitors of *sn*-1-diacylglycerol lipase  $\alpha$  and  $\alpha/\beta$ -hydrolase domain 6. *J. Med. Chem.* **57**(15), 6610-6622 (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, 02/09/2024

#### 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