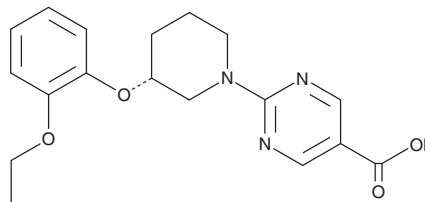


# PRODUCT INFORMATION

## (R)-2-(3-(2-Ethoxyphenoxy)piperidin-1-yl)pyrimidine-5-carboxylic Acid

Item No. 39257

CAS Registry No.: 1809064-89-8  
 Formal Name: 2-[(3R)-3-(2-ethoxyphenoxy)-1-piperidinyl]-5-pyrimidinecarboxylic acid  
 MF:  $C_{18}H_{21}N_3O_4$   
 FW: 343.4  
 Purity:  $\geq 95\%$   
 Supplied as: A solid  
 Storage:  $-20^{\circ}\text{C}$   
 Stability:  $\geq 4$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Description

(R)-2-(3-(2-Ethoxyphenoxy)piperidin-1-yl)pyrimidine-5-carboxylic acid is an intermediate in the synthesis of the diacylglycerol acyltransferase 2 (DGAT-2) inhibitor ervogastat (Item No. 37929).<sup>1</sup>

### Reference

1. Futatsugi, K., Cabral, S., Kung, D.W., *et al.* Discovery of ervogastat (PF-06865571): A potent and selective inhibitor of diacylglycerol acyltransferase 2 for the treatment of non-alcoholic steatohepatitis. *J. Med. Chem.* **65**(22), 15000-15013 (2022).

#### 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, 08/11/2023

#### 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