

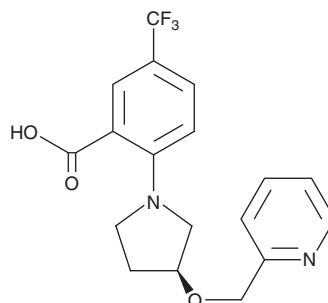
# PRODUCT INFORMATION



## XEN445

Item No. 19144

**CAS Registry No.:** 1515856-92-4  
**Formal Name:** 2-[(3S)-3-(2-pyridinylmethoxy)-1-pyrrolidinyl]-5-(trifluoromethyl)-benzoic acid  
**MF:** C<sub>18</sub>H<sub>17</sub>F<sub>3</sub>N<sub>2</sub>O<sub>3</sub>  
**FW:** 366.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 274, 329 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

XEN445 is supplied as a crystalline solid. A stock solution may be made by dissolving the XEN445 in the solvent of choice, which should be purged with an inert gas. XEN445 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of XEN445 in these solvents is approximately 14, 25, and 30 mg/ml, respectively.

XEN445 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, XEN445 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. XEN445 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

Endothelial lipase, a member of the triglyceride lipase gene family expressed in endothelial cells, prefers phospholipid-enriched high-density lipoprotein (HDL) over triglyceride as a substrate. XEN445 is an endothelial lipase inhibitor (IC<sub>50</sub> = 237 nM) that demonstrates selectivity for endothelial lipase compared to lipoprotein and hepatic lipases (IC<sub>50</sub>s = 20 and 9.5 μM, respectively).<sup>1</sup> An oral dose of 30 mg/kg XEN445 is reported to increase HDL cholesterol concentrations by 16% after three days and by 30% after nine days dosing in wild-type mice.<sup>1</sup>

### Reference

1. Sun, S., Dean, R., Jia, Q., *et al.* Discovery of XEN445: A potent and selective endothelial lipase inhibitor raises plasma HDL-cholesterol concentration in mice. *Bioorg. Med. Chem.* **21(24)**, 7724-7734 (2013).

#### 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, 11/22/2022

#### 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](http://WWW.CAYMANCHEM.COM)