

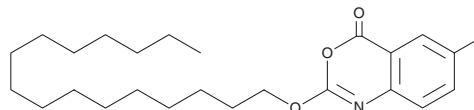
# PRODUCT INFORMATION



## Cetilistat

Item No. 31622

**CAS Registry No.:** 282526-98-1  
**Formal Name:** 2-(hexadecyloxy)-6-methyl-4H-3,1-benzoxazin-4-one  
**Synonym:** ATL-962  
**MF:** C<sub>25</sub>H<sub>39</sub>NO<sub>3</sub>  
**FW:** 401.6  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 223, 228 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

Cetilistat is supplied as a solid. A stock solution may be made by dissolving the cetilistat in the solvent of choice, which should be purged with an inert gas. Cetilistat is slightly soluble in acetonitrile and DMSO.

### Description

Cetilistat is an inhibitor of pancreatic lipase (IC<sub>50</sub>s = 5.95 and 54.8 nM for the human and rat enzymes, respectively).<sup>1</sup> It reduces increases in plasma triglyceride levels in an oral fat tolerance test, indicating decreased intestinal fat absorption, in rats when administered at a doses ranging from 3 to 100 mg/kg. Dietary administration of cetilistat (4.9-50.7 mg/kg per day) decreases body weight gain without affecting food intake, as well as reduces increases in plasma triglyceride and leptin levels in a rat model of high-fat diet-induced obesity. It also reduces viral load in the culture supernatant of Vero E6 cells infected with severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2; EC<sub>90</sub> = 2.9 μM) and inhibits SARS-CoV-2 replication in a plaque reduction assay (EC<sub>50</sub> = 1.13 μM).<sup>2</sup>

### References

1. Yamada, Y., Kato, T., Ogino, H., *et al.* Cetilistat (ATL-962), a novel pancreatic lipase inhibitor, ameliorates body weight gain and improves lipid profiles in rats. *Horm. Metab. Res.* **40(8)**, 539-543 (2008).
2. Yuan, S., Chan, J.F.W., Chik, K.K.H., *et al.* Discovery of the FDA-approved drugs bexarotene, cetilistat, diiodohydroxyquinoline, and abiraterone as potential COVID-19 treatments with a robust two-tier screening system. *Pharmacol. Res.* **159**, 104960 (2020).

#### 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, 10/28/2022

#### 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