

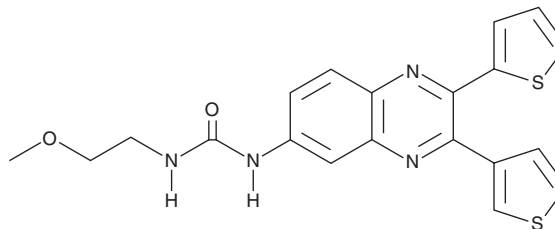
# PRODUCT INFORMATION



## Acetyl-CoA Synthase Inhibitor

Item No. 36023

**CAS Registry No.:** 508186-14-9  
**Formal Name:** N-(2,3-di-2-thienyl-6-quinoxaliny)-N'-(2-methoxyethyl)-urea  
**Synonyms:** ACCS2 Inhibitor, MDK6149, VY-3-249  
**MF:** C<sub>20</sub>H<sub>18</sub>N<sub>4</sub>O<sub>2</sub>S<sub>2</sub>  
**FW:** 410.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 221, 273, 400 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

Acetyl-CoA synthase inhibitor is supplied as a solid. A stock solution may be made by dissolving the acetyl-CoA synthase inhibitor in the solvent of choice, which should be purged with an inert gas. Acetyl-CoA synthase inhibitor is soluble in DMSO.

### Description

Acetyl-CoA synthase inhibitor is an inhibitor of acetyl-CoA synthetase 2 (ACCS2; IC<sub>50</sub> = 1.214 μM).<sup>1</sup> It reduces oleic acid-induced increases in the number of histone H2AX-positive cells, a marker of DNA damage, in cultured human immortalized hepatocytes when used at a concentration of 50 μM.<sup>2</sup>

### References

- Schug, Z. and Salvino, J. Acetyl-CoA synthetase 2 (ACCS2) inhibitors and methods using same. *The Wistar Institute*. **WO2020/252407A1** (2020).
- Assante, G., Chandrasekaran, S., Ng, S., *et al.* Acetyl-CoA metabolism drives epigenome change and contributes to carcinogenesis risk in fatty liver disease. *Genome Med.* **14(1)**, 67 (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, 12/19/2022

#### 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)