

# PRODUCT INFORMATION



**NB-598**

Item No. 14912

**CAS Registry No.:** 131060-14-5

**Formal Name:** 3-([3,3'-bithiophen]-5-ylmethoxy)-N-[(2E)-6,6-dimethyl-2-hepten-4-yn-1-yl]-N-ethyl-benzenemethanamine

**MF:** C<sub>27</sub>H<sub>31</sub>NOS<sub>2</sub>

**FW:** 449.7

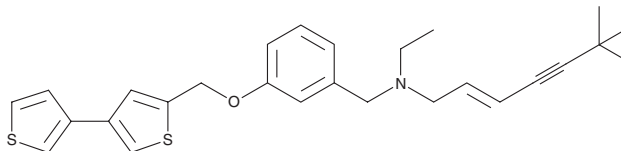
**Purity:** ≥98%

**UV/Vis.:** λ<sub>max</sub>: 219 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

NB-598 is supplied as a solid. A stock solution may be made by dissolving the NB-598 in the solvent of choice, which should be purged with an inert gas. NB-598 is soluble in ethanol (heated) and DMSO (heated).

## Description

NB-598 is an inhibitor of squalene epoxidase (IC<sub>50</sub>s = 2 and 0.75 nM for dog liver and human microsomal enzymes, respectively).<sup>1</sup> It selectively inhibits cholesterol synthesis over free fatty acid, triglyceride, and phospholipid synthesis in HepG2 cells at 1 μM. NB-598 (10 mg/kg per day) reduces serum cholesterol levels in dogs.

## Reference

1. Horie, M., Tsuchiya, Y., Hayashi, M., *et al.* NB-598: A potent competitive inhibitor of squalene epoxidase. *J. Biol. Chem.* **265**(30), 18075-18078 (1990).

### 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, 05/02/2023

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