

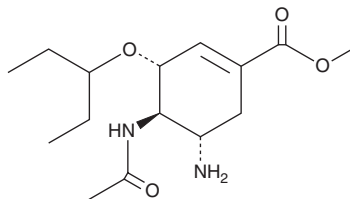
# PRODUCT INFORMATION



## Oseltamivir Acid methyl ester

Item No. 36556

**CAS Registry No.:** 208720-71-2  
**Formal Name:** (3R,4R,5S)-4-(acetylamino)-5-amino-3-(1-ethylpropoxy)-1-cyclohexene-1-carboxylic acid, methyl ester  
**Synonym:** ZINC036451498  
**MF:**  $C_{15}H_{26}N_2O_4$   
**FW:** 298.4  
**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.

### Laboratory Procedures

Oseltamivir acid methyl ester is supplied as a solid. A stock solution may be made by dissolving the oseltamivir acid methyl ester in the solvent of choice, which should be purged with an inert gas. Oseltamivir acid methyl ester is slightly soluble in chloroform and methanol.

### Description

Oseltamivir acid methyl ester is a prodrug form of the neuraminidase inhibitor and antiviral oseltamivir acid (Item No. 15779).<sup>1</sup> It is converted to oseltamivir acid by carboxylesterase 1 (CES1).

### Reference

1. Shimizu, M., Fukami, T., Ogawa, H., *et al.* Systematic approach for screening of prodrugs: Evaluation using oseltamivir analogues as models. *J. Pharm. Sci.* **110**(2), 925-934 (2021).

#### 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, 12/20/2022

#### 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)