

PRODUCT INFORMATION



Laninamivir octanoate

Item No. 40850

CAS Registry No.: 203120-46-1

Formal Name: 5-(acetylamino)-4-
[(aminoiminomethyl)amino]-2,6-
anhydro-3,4,5-trideoxy-7-O-
methyl-D-glycero-D-galacto-non-
2-enonic acid, 9-octanoate

Synonym: CS 8958

MF: $C_{21}H_{36}N_4O_8$

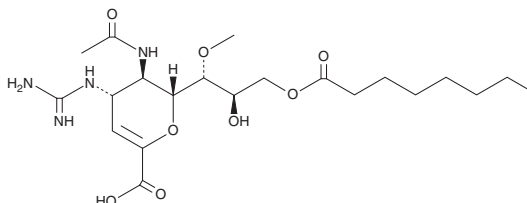
FW: 472.5

Purity: $\geq 98\%$

Supplied as: A solid

Storage: -20°C

Stability: ≥ 2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Laninamivir octanoate is supplied as a solid. A stock solution may be made by dissolving the laninamivir octanoate in the solvent of choice, which should be purged with an inert gas. Laninamivir octanoate is slightly soluble (0.1-1 mg/ml) in ethanol and sparingly soluble (1-10 mg/ml) in DMSO.

Description

Laninamivir octanoate is an inhibitor of influenza neuraminidases and a prodrug form of laninamivir.¹ It selectively inhibits the neuraminidases from influenza A H1N1, H2N2, and H3N2 strains (mean IC_{50} s = 822, 128, and 114 nM, respectively) over the neuraminidases from influenza B Victoria and Yamagata strains (mean IC_{50} s = $>10,000$ nM for both). Intranasal administration of laninamivir octanoate (0.95 $\mu\text{g/kg}$) increases survival in a mouse model of lethal influenza H1N1 infection. Formulations containing laninamivir octanoate have been used in the prevention of influenza.

Reference

1. Yamashita, M., Tomozawa, T., Kakuta, M., *et al.* CS-8958, a prodrug of the new neuraminidase inhibitor R-125489, shows long-acting anti-influenza virus activity. *Antimicrob. Agents Chemother.* **53**(1), 186-192 (2009).

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, 04/15/2024

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