

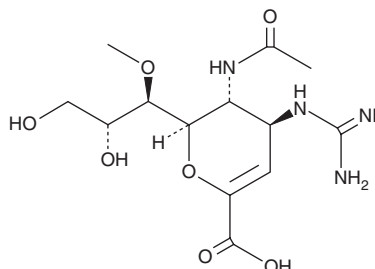
PRODUCT INFORMATION



Laninamivir

Item No. 39139

CAS Registry No.: 203120-17-6
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
Synonym: R-125489
MF: C₁₃H₂₂N₄O₇
FW: 346.3
Purity: ≥98%
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

Laninamivir is supplied as a solid. Aqueous solutions of laninamivir can be prepared by directly dissolving the solid in aqueous buffers. Laninamivir is sparingly soluble (1-10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Laninamivir is an inhibitor of influenza neuraminidases and an active metabolite of the prodrug laninamivir octanoate (Item No. 40850).¹ It inhibits the neuraminidases from influenza A H1N1, H2N2, and H3N2 strains (mean IC₅₀s = 2.49, 11.4, and 16.7 nM, respectively) and influenza B Victoria and Yamagata strains (mean IC₅₀s = 18.9 nM for both). Intranasal administration of laninamivir (69 µg/kg) increases survival in a mouse model of lethal influenza H1N1 infection. Formulations containing laninamivir 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, 08/02/2024

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