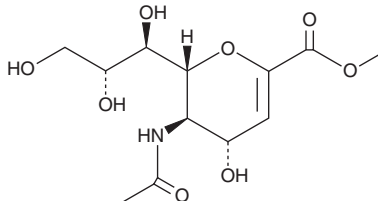


PRODUCT INFORMATION

N-acetyl-2,3-dehydro-2-Deoxyneuraminic Acid methyl ester

Item No. 36529

CAS Registry No.: 25875-99-4
Formal Name: 5-(acetylamino)-2,6-anhydro-3,5-dideoxy-D-glycero-D-galacto-non-2-enonic acid, methyl ester
MF: C₁₂H₁₉NO₈
FW: 305.3
Purity: ≥95%
UV/Vis.: λ_{max}: 245 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

N-acetyl-2,3-dehydro-2-Deoxyneuraminic acid methyl ester is supplied as a solid. Aqueous solutions of N-acetyl-2,3-dehydro-2-deoxyneuraminic acid methyl ester can be prepared by directly dissolving the solid in aqueous buffers. The solubility of N-acetyl-2,3-dehydro-2-deoxyneuraminic acid methyl ester in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

N-acetyl-2,3-dehydro-2-Deoxyneuraminic acid methyl ester is a bacterial neuraminidase inhibitor (K_i = 48 μM for the *A. sialophilus* enzyme).¹

Reference

- Flashner, M., Kessler, J., and Tanenbaum, S.W. The interaction of N-acetylneuraminic acid and its analogs with neuraminidase. *Develop. Biochem.* 6, 27-44 (1979).

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, 05/05/2023

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