

PRODUCT INFORMATION



Peramivir

Item No. 23765

CAS Registry No.: 330600-85-6
Formal Name: (1S,2S,3R,4R)-3-[(1S)-1-(acetylamino)-2-ethylbutyl]-4-[(aminoiminomethyl)amino]-2-hydroxy-cyclopentanecarboxylic acid

Synonyms: BCX 1812, RWJ 270201

MF: C₁₅H₂₈N₄O₄

FW: 328.4

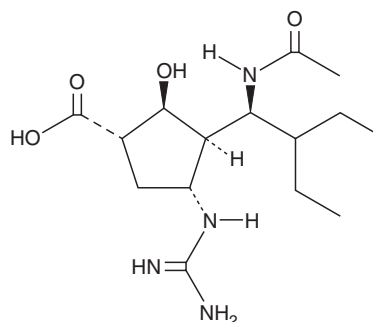
Purity: ≥98%

Supplied as: A crystalline 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

Peramivir is supplied as a crystalline solid. A stock solution may be made by dissolving the peramivir in the solvent of choice, which should be purged with an inert gas. Peramivir is slightly soluble in ethanol and methanol.

Description

Peramivir is an inhibitor of influenza neuraminidase (IC₅₀s = 0.09 and 11 nM for influenza A and B neuraminidases, respectively).¹ It is selective for influenza neuraminidase over bacterial, mammalian, and other viral neuraminidases (IC₅₀s = >300 μM). Peramivir inhibits neuraminidase activity in H1N1, H2N2, H3N2, and H6N2 influenza strains (IC₅₀s = 0.09-1.1 nM) and reduces lysis of MDCK cells infected with influenza (EC₅₀s = <0.01-21 nM). Pretreatment with peramivir (10-100 mg/kg) protects mice against lethal influenza infections. It also increases survival in ferrets infected intranasally with avian influenza type A H5N1 when injected intramuscularly after infection.² Formulations containing peramivir have been used to treat influenza.

References

1. Young, D., Fowler, C., and Bush, K. RWJ-270201 (BCX-1812): A novel neuraminidase inhibitor for influenza. *Philos. Trans. R. Soc. Lond. B Biol. Sci* **356(1416)**, 1905-1913 (2001).
2. Yun, N.E., Linde, N.S., Zacks, M.A., et al. Injectable peramivir mitigates disease and promotes survival in ferrets and mice infected with the highly virulent influenza virus, A/Vietnam/1203/04 (H5N1). *Virology* **374(1)**, 198-209 (2008).

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, 12/05/2022

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