

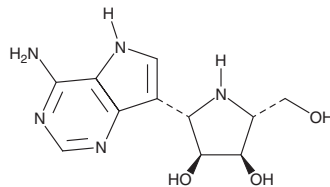
PRODUCT INFORMATION



BCX4430

Item No. 16085

CAS Registry No.: 249503-25-1
Formal Name: (2S,3S,4R,5R)-2-(4-amino-5H-pyrrolo[3,2-d]pyrimidin-7-yl)-5-(hydroxymethyl)-3,4-pyrrolidinediol
Synonym: Galidesivir
MF: C₁₁H₁₅N₅O₃
FW: 265.3
Purity: ≥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

BCX4430 is supplied as a solid. A stock solution may be made by dissolving the BCX4430 in water. The solubility of BCX4430 in water is approximately 1.5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

BCX4430 is an adenosine nucleoside analog with broad-spectrum antiviral activity with EC₅₀ values ranging from 3.4 to >100 μM against a panel of Ebola, Marburg, bunya-, arena-, paramyxo-, corona- and flaviviruses.¹ It undergoes anabolism to BCX4430-triphosphate in Huh-7 cells followed by pyrophosphate cleavage to BCX4430-monophosphate, which is incorporated into RNA strands to induce non-obligate RNA chain termination. BCX4430 inhibits Ebola virus minigenome RNA replication and hepatitis C virus (HCV) NS5B RNA polymerase activity in a dose-dependent manner. *In vivo*, BCX4430 (5-150 mg/kg) increases survival in mice infected with Rift Valley fever virus. It increases survival of Marburg virus-Musoke infected guinea pigs and cynomolgus macaques when administered pre- or post-infection. BCX4430 (300 mg/kg) also protects mice from mortality in a mouse model of severe Zika virus infection.²

References

1. Warren, T.K., Wells, J., Panchal, R.G., *et al.* Protection against filovirus diseases by a novel broad-spectrum nucleoside analogue BCX4430. *Nature* **508**(7496), 402-405 (2014).
2. Julander, J.G., Siddharthan, V., Evans, J., *et al.* Efficacy of the broad-spectrum antiviral compound BCX4430 against Zika virus in cell culture and in a mouse model. *Antiviral Res.* **137**, 14-22 (2017).

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, 03/24/2020

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