

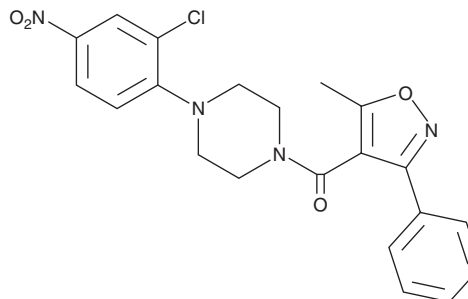
PRODUCT INFORMATION



Nucleozin

Item No. 30363

CAS Registry No.: 341001-38-5
Formal Name: [4-(2-chloro-4-nitrophenyl)-1-piperazinyl](5-methyl-3-phenyl-4-isoxazolyl)-methanone
MF: C₂₁H₁₉ClN₄O₄
FW: 426.9
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

Nucleozin is supplied as a crystalline solid. A stock solution may be made by dissolving the nucleozin in the solvent of choice, which should be purged with an inert gas. Nucleozin is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of nucleozin in these solvents is approximately 30 mg/ml.

Description

Nucleozin is an antiviral inhibitor of influenza nucleoprotein accumulation, preventing it from translocating into the nucleus of host cells.¹ It inhibits MDCK cell infection by influenza A and H5N1 (EC₅₀s = 0.069 and 0.33 μM, respectively), as well as a clinical isolate of H3N2 (EC₅₀ = 0.16 μM). Nucleozin reduces the lung viral load and increases the survival rate in a mouse model of influenza H5N1 infection.

Reference

1. Kao, R.Y., Yang, D., Lau, L.-S., *et al.* Identification of influenza A nucleoprotein as an antiviral target. *Nat. Biotechnol.* **26**(6), 600-605 (2010).

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