

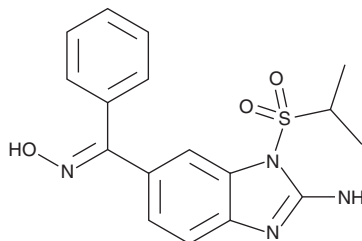
PRODUCT INFORMATION



Enviroxime

Item No. 30552

CAS Registry No.: 72301-79-2
Formal Name: (1E)-[2-amino-1-[(1-methylethyl)sulfonyl]-1H-benzimidazol-6-yl]phenyl-methanone oxime
Synonym: LY122772
MF: C₁₇H₁₈N₄O₃S
FW: 358.4
Purity: ≥95%
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

Enviroxime is supplied as a solid. A stock solution may be made by dissolving the enviroxime in the solvent of choice, which should be purged with an inert gas. Enviroxime is soluble in DMSO. Enviroxime is slightly soluble in methanol (warmed).

Description

Enviroxime is an antiviral agent.¹ It inhibits coxsackievirus B3-, rhinovirus-, or poliovirus-induced plaque formation in HeLa monolayer cell cultures when used at concentrations of 0.1, 1, and 10 µg/ml.¹ Enviroxime (1 µg/ml) inhibits production of poliovirus PV1 viral RNA in HeLa monolayer cell cultures. It reduces viral titers in DRAW macrophages infected with the DA strain of Theiler's murine encephalomyelitis virus.²

References

1. Heinz, B.A. and Vance, L.M. The antiviral compound enviroxime targets the 3A coding region of rhinovirus and poliovirus. *J. Virol.* **69**(7), 4189-4197 (1995).
2. Steurbaut, S., Merckx, E., Rombaut, B., *et al.* Modulation of viral replication in macrophages persistently infected with the DA strain of Theiler's murine encephalomyelitis virus. *Virol. J.* **5**, 89 (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, 05/19/2020

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