

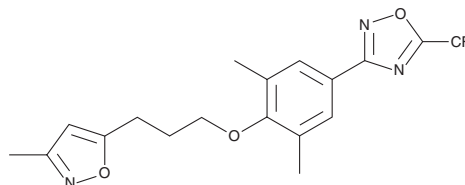
# PRODUCT INFORMATION



## Pleconaril

Item No. 28461

**CAS Registry No.:** 153168-05-9  
**Formal Name:** 3-[3,5-dimethyl-4-[3-(3-methyl-5-isoxazolyl)propoxy]phenyl]-5-(trifluoromethyl)-1,2,4-oxadiazole  
**Synonyms:** VP 63843, WIN 63,843  
**MF:** C<sub>18</sub>H<sub>18</sub>F<sub>3</sub>N<sub>3</sub>O<sub>3</sub>  
**FW:** 381.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 262 nm  
**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

Pleconaril is supplied as a crystalline solid. A stock solution may be made by dissolving the pleconaril in the solvent of choice, which should be purged with an inert gas. Pleconaril is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of pleconaril in these solvents is approximately 10, 20, and 25 mg/ml, respectively.

### Description

Pleconaril is an antipicornaviral agent.<sup>1,2</sup> It inhibits *Picornaviridae* viral replication by binding to a hydrophobic pocket in the major VP1 capsid protein, which prevents uncoating of the viral RNA genome. Pleconaril inhibits replication of the rhinoviruses HRV-A2 and HRV-B14 in HeLa Rh cells (EC<sub>50</sub>s = 0.1 and 0.3 μM, respectively) and is not cytotoxic to HeLa Rh cells with a 50% cytotoxic concentration (CC<sub>50</sub>) of greater than 131 μM.<sup>2</sup> It also inhibits replication of enterovirus 71 (EV71) in human rhabdomyosarcoma (RD) cells (EC<sub>50</sub> = 15 μM) and is not cytotoxic to RD cells (CC<sub>50</sub> = >131 μM). Pleconaril (80 mg/kg per day) increases survival of EV71-infected mouse pups from 20 to 80%.<sup>3</sup>

### References

1. Florea, N.R., Maglio, D., and Nicolau, D.P. Pleconaril, a novel antipicornaviral agent. *Pharmacotherapy* **23**(3), 339-348 (2003).
2. Bernard, A., Lacroix, C., Cabiddu, M.G., et al. Exploration of the anti-enterovirus activity of a series of pleconaril/pirodavir-like compounds. *Antivir. Chem. Chemother.* **24**(2), 56-61 (2015).
3. Zhang, G., Zhou, F., Gu, B., et al. *In vitro* and *in vivo* evaluation of ribavirin and pleconaril antiviral activity against enterovirus 71 infection. *Arch. Virol.* **157**(4), 669-679 (2012).

#### 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, 06/26/2024

#### 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](http://WWW.CAYMANCHEM.COM)