

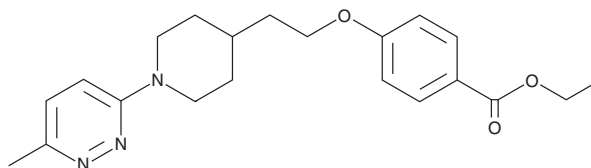
# PRODUCT INFORMATION



## Pirodavir

Item No. 22922

**CAS Registry No.:** 124436-59-5  
**Formal Name:** 4-[2-[1-(6-methyl-3-pyridazinyl)-4-piperidinyl]ethoxy]benzoic acid, ethyl ester  
**Synonym:** R-77975  
**MF:** C<sub>21</sub>H<sub>27</sub>N<sub>3</sub>O<sub>3</sub>  
**FW:** 369.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 257 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

Pirodavir is supplied as a crystalline solid. A stock solution may be made by dissolving the pirodavir in the solvent of choice, which should be purged with an inert gas. Pirodavir is soluble in the organic solvent dimethyl formamide (DMF) at a concentration of approximately 30 mg/ml.

Pirodavir is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, pirodavir should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Pirodavir has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Pirodavir is a broad-spectrum antipicornaviral agent.<sup>1</sup> It is active against 80 typed (MICs = 0.001-14.4 µg/ml) and one untyped rhinovirus strain (MIC = 0.002 µg/ml) and is active against both group A and group B serotypes. It is also active against serotypes of coxsackievirus, poliovirus, and echovirus enteroviruses (MICs = 0.018-8.5, 0.025-0.27, and 0.019-1.3 µg/ml, respectively).

### Reference

1. Andries, K., Dewindt, B., Snoeks, J., *et al.* In vitro activity of pirodavir (R 77975), a substituted phenoxy-pyridazinamine with broad-spectrum antipicornaviral activity. *Antimicrob. Agents Chemother.* **36**(1), 100-107 (1992).

#### 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, 11/04/2022

#### 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