

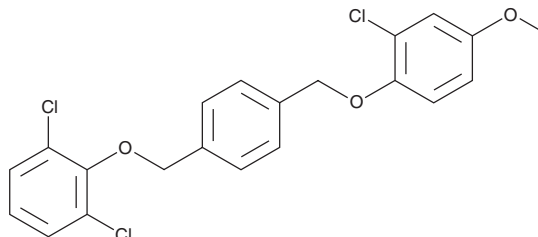
# PRODUCT INFORMATION



## Pocapavir

Item No. 37349

**CAS Registry No.:** 146949-21-5  
**Formal Name:** 1-[(2-chloro-4-methoxyphenoxy)methyl]-4-[(2,6-dichlorophenoxy)methyl]-benzene  
**Synonyms:** SCH 48973, V-073  
**MF:** C<sub>21</sub>H<sub>17</sub>Cl<sub>3</sub>O<sub>3</sub>  
**FW:** 423.7  
**Purity:** ≥98%  
**Supplied as:** A 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

Pocapavir is supplied as a solid. A stock solution may be made by dissolving the pocapavir in the solvent of choice, which should be purged with an inert gas. Pocapavir is soluble in organic solvents such as methanol and DMSO.

### Description

Pocapavir is an antiviral agent.<sup>1</sup> It reduces viral replication in HeLa cells infected with poliovirus type 2, group A9 coxsackievirus, or Echovirus type 4, -6, or -11 (IC<sub>50</sub>s = 0.08, 0.09, 0.11, 0.03, and 0.02 µg/ml, respectively). It also inhibits viral replication of 154 clinical enterovirus isolates in infected HeLa cells (mean IC<sub>50</sub> = 0.9 µg/ml). *In vivo*, pocapavir (3, 10, and 20 mg/kg) increases survival in poliovirus type 2-infected mice.

### Reference

1. Buontempo, P.J., Cox, S., Wright-Minogue, J., *et al.* SCH 48973: A potent, broad-spectrum, antienterovirus compound. *Antimicrob Agents Chemother.* **41(6)**, 1220-1225 (1997).

#### 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, 11/01/2023

#### 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