

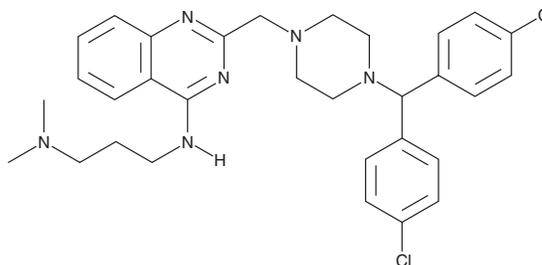
# PRODUCT INFORMATION



**SCH 529074**

Item No. 21552

**CAS Registry No.:** 922150-11-6  
**Formal Name:** N<sup>3</sup>-[2-[[4-[bis(4-chlorophenyl)methyl]-1-piperazinyl]methyl]-4-quinazoliny]-N<sup>1</sup>,N<sup>1</sup>-dimethyl-1,3-propanediamine  
**MF:** C<sub>31</sub>H<sub>36</sub>Cl<sub>2</sub>N<sub>6</sub>  
**FW:** 563.6  
**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

SCH 529074 is supplied as a crystalline solid. A stock solution may be made by dissolving the SCH 529074 in the solvent of choice. SCH 529074 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of SCH 529074 in DMSO is approximately 25 mg/ml and approximately 33 mg/ml in ethanol and DMF.

## Description

SCH 529074 is a small molecule activator of p53, a tumor suppressor that is mutated or non-functional in 50% of human tumors.<sup>1</sup> SCH 529074 restores DNA binding activity and increases DNA binding affinity of two recombinant GST fusion mutant p53 DNA binding domains (R273H and R249S) in an electrophoretic mobility shift assay with K<sub>d</sub> values of 0.6740 and 0.4313, respectively. It also stimulates DNA binding activity and prevents human double minute-2 (Hdm-2) ubiquitination of wild-type p53 *in vitro*. SCH 29074 reduces proliferation of tumor cell lines with mutant and wild-type p53 (EC<sub>50</sub>s = 400-3,700 nM), with the p53 mutant cell lines being the most sensitive. *In vivo*, SCH 529074 reduces tumor growth in a DLD-1 human colorectal cancer xenograft model in mice in a dose-dependent manner.

## Reference

1. Demma, M., Maxwell, E., Ramos, R., *et al.* SCH529074, a small molecule activator of mutant p53, which binds p53 DNA binding domain (DBD), restores growth-suppressive function to mutant p53 and interrupts HDM2-mediated ubiquitination of wild type p53. *J. Biol. Chem.* **285**(14), 10198-10212 (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, 12/22/2022

## 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