

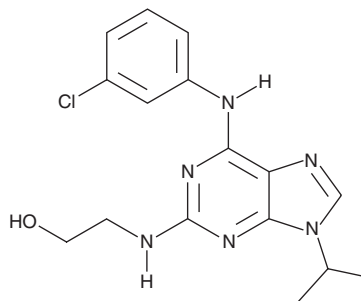
PRODUCT INFORMATION



NG 52

Item No. 17309

CAS Registry No.: 212779-48-1
Formal Name: 2-[[6-[(3-chlorophenyl)amino]-9-(1-methylethyl)-9H-purin-2-yl]amino]-ethanol
Synonym: Compound 52
MF: C₁₆H₁₉ClN₆O
FW: 346.8
Purity: ≥98%
UV/Vis.: λ_{max}: 215, 260, 313 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

NG 52 is supplied as a crystalline solid. A stock solution may be made by dissolving the NG 52 in the solvent of choice, which should be purged with an inert gas. NG 52 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of NG 52 in ethanol is approximately 5 mg/ml and approximately 15 mg/ml in DMSO and DMF.

NG 52 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, NG 52 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. NG 52 has a solubility of approximately 0.1 mg/ml in a 1:7 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

NG 52 is a tri-substituted purine that binds to the ATP-binding site of yeast cyclin-dependent kinases, inhibiting Cdc28p and Pho85p (IC₅₀s = 7 and 2 μM, respectively).¹ It is ineffective against the yeast kinases Kin28p, Srb10, and Cak1p. NG 52 is cell permeable and inhibits the growth of *S. cerevisiae* (GI₅₀ = 30 μM).¹ It is an analog of purvalanol A (Item No. 14579), a potent inhibitor of mammalian cyclin-dependent kinases.¹

Reference

1. Gray, N.S., Wodicka, L., Thunnissen, A.M.W.H., et al. Exploiting chemical libraries, structure, and genomics in the search for kinase inhibitors. *Science* 281(5376), 533-538 (1998).

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, 10/03/2022

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