

PRODUCT INFORMATION



RO8994

Item No. 25674

CAS Registry No.: 1309684-94-3
Formal Name: (2'S,3R,4'S,5'R)-N-[4-(aminocarbonyl)-2-methoxyphenyl]-6-chloro-4'-(3-chloro-2-fluorophenyl)-2'-(2,2-dimethylpropyl)-1,2-dihydro-2-oxo-spiro[3H-indole-3,3'-pyrrolidine]-5'-carboxamide

MF: C₃₁H₃₁Cl₂FN₄O₄

FW: 613.5

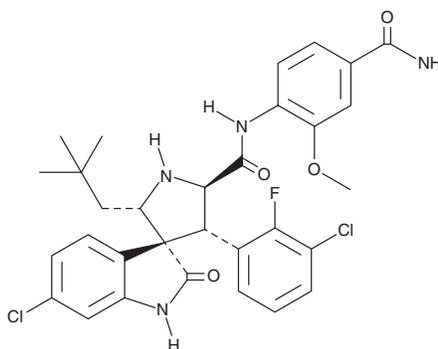
Purity: ≥98%

UV/Vis.: λ_{max}: 212, 270, 300 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

RO8994 is supplied as a crystalline solid. A stock solution may be made by dissolving the RO8994 in the solvent of choice, which should be purged with an inert gas. RO8994 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of RO8994 in these solvents is approximately 5, 14, and 12.5 mg/ml, respectively.

Description

RO8994 is an inhibitor of the protein-protein interaction between MDM2 and the tumor suppressor p53 in a homogeneous time-resolved fluorescence (HTRF) binding assay (IC₅₀ = 5 nM).¹ It inhibits the proliferation of SJSA-1, RKO, and HCT116 cells with an average IC₅₀ value of 20 nM and increases apoptosis in SJSA-1 cells in a concentration-dependent manner. RO8994 reduces tumor growth and induces tumor regression in an SJSA-1 osteosarcoma mouse xenograft model when administered at doses of 1.56 and 6.25 mg/kg, respectively.

Reference

1. Zhang, Z., Ding, Q., Liu, J.J., *et al.* Discovery of potent and selective spiroindolinone MDM2 inhibitor, RO8994, for cancer therapy. *Bioorg. Med. Chem.* **22(15)**, 4001-4009 (2014).

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, 11/28/2022

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