

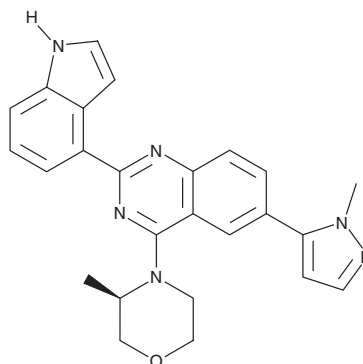
# PRODUCT INFORMATION



## SKLB197

Item No. 45728

**CAS Registry No.:** 2713577-16-1  
**Formal Name:** 2-(1H-indol-4-yl)-4-[(3R)-3-methyl-4-morpholinyl]-6-(1methyl-1H-pyrazol-5-yl)-quinazoline  
**MF:** C<sub>25</sub>H<sub>24</sub>N<sub>6</sub>O  
**FW:** 424.5  
**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

SKLB197 is supplied as a solid. A stock solution may be made by dissolving the SKLB197 in the solvent of choice, which should be purged with an inert gas. SKLB197 is sparingly soluble (1-10 mg/ml) in DMSO.

### Description

SKLB197 is an inhibitor of ataxia-telangiectasia and Rad3-related protein/kinase (ATR) ( $IC_{50} = 0.013 \mu M$ ).<sup>1</sup> It is selective for ATR over a panel of 402 other kinases ( $IC_{50}s = >1 \mu M$ ). SKLB197 reduces growth in cancer cell lines deficient in ataxia-telangiectasia mutated kinase (ATM;  $IC_{50}s = 0.13-0.66 \mu M$ ) to a greater extent than in ATM-expressing cancer cell lines ( $IC_{50}s = 3.66-8.79 \mu M$ ). It induces apoptosis in ATM-deficient LoVo colorectal cancer cells but not in ATM-expressing HT-29 colorectal cancer cells and reduces tumor growth in a LoVo mouse xenograft model when administered at doses of 50 or 100 mg/kg.

### Reference

1. Bin, H., Chen, P., Wu, M., *et al.* Discovery of a potent and highly selective inhibitor of ataxia telangiectasia mutated and Rad3-Related (ATR) kinase: Structural activity relationship and antitumor activity both *in vitro* and *in vivo*. *Eur. J. Med. Chem.* **232**, 114187 (2022).

#### 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, 05/26/2026

#### 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