

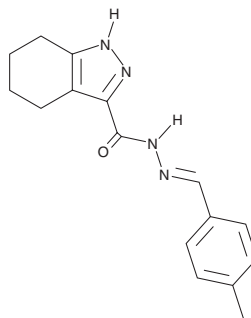
# PRODUCT INFORMATION



## Suprafenacine

Item No. 31158

**CAS Registry No.:** 1477482-50-0  
**Formal Name:** 4,5,6,7-tetrahydro-1H-indazole-3-carboxylic acid (2E)-2-[(4-methylphenyl)methylene]hydrazide  
**Synonym:** SML1493  
**MF:** C<sub>16</sub>H<sub>18</sub>N<sub>4</sub>O  
**FW:** 282.3  
**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

Suprafenacine is supplied as a solid. A stock solution may be made by dissolving the suprafenacine in the solvent of choice, which should be purged with an inert gas. Suprafenacine is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml.

### Description

Suprafenacine is an inhibitor of tubulin polymerization ( $IC_{50} = 0.38 \mu M$ ) with anticancer activity.<sup>1</sup> It inhibits the proliferation of HeLa, Jurkat, SNU-16, MDA-MB-231, A549, SH-SY5Y, HCT15, and KB-3-1 cells ( $IC_{50}$ s = 83-381.2 nM) and P-glycoprotein-enriched KB-VI cells ( $IC_{50} = 215.7$  nM). Suprafenacine (10  $\mu M$ ) increases phosphorylation of JNK and Bcl-2 and induces apoptosis and cell cycle arrest at the G<sub>2</sub>/M phase in HeLa cells. Suprafenacine (20 mg/kg) reduces tumor growth in an HCT15 colon adenocarcinoma mouse xenograft model.

### Reference

1. Choi, B.-H., Chattopadhyaya, S., Thanh, L.N., *et al.* Suprafenacine, an indazole-hydrazide agent, targets cancer cells through microtubule destabilization. *PLoS One* **9(10)**, e110955 (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, 11/03/2022

#### 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](http://WWW.CAYMANCHEM.COM)