

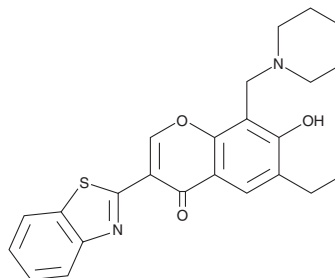
# PRODUCT INFORMATION



## SZL P1-41

Item No. 26142

**CAS Registry No.:** 222716-34-9  
**Formal Name:** 3-(2-benzothiazolyl)-6-ethyl-7-hydroxy-8-(1-piperidinylmethyl)-4H-1-benzopyran-4-one  
**MF:** C<sub>24</sub>H<sub>24</sub>N<sub>2</sub>O<sub>3</sub>S  
**FW:** 420.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 224, 264, 308 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

SZL P1-41 is supplied as a crystalline solid. A stock solution may be made by dissolving the SZL P1-41 in the solvent of choice, which should be purged with an inert gas. SZL P1-41 is soluble in ethanol, DMSO, and dimethyl formamide (DMF). SZL P1-41 has a solubility of approximately 0.3 mg/ml in ethanol and DMSO and approximately 1 mg/ml in DMF.

### Description

SZL P1-41 is an inhibitor of the Skp1-Cullin1-F-box (SCF) family protein Skp2.<sup>1</sup> It inhibits the interaction between Skp2 and Skp1 to block Skp2 E3 ubiquitin ligase activity when used at a concentration of 5 μM. SZL P1-41 selectively inhibits Skp2 over Fbw7 and β-TrCP E3 ubiquitin ligase activity in HEK293T cells. It reduces survival of LNCaP and PC3 prostate cancer cells in a concentration-dependent manner via inhibition of aerobic glycolysis and induction of apoptosis. SZL P1-41 (40 and 80 mg/kg) reduces tumor volume in A549 lung cancer and PC3 prostate cancer mouse xenograft models.

### Reference

1. Chan, C.H., Morrow, J.K., Li, C.F., *et al.* Pharmacological inactivation of Skp2 SCF ubiquitin ligase restricts cancer stem cell traits and cancer progression. *Cell*. **154**(3), 556-568 (2013).

#### 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, 12/01/2022

#### 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