

PRODUCT INFORMATION



ML-60218

Item No. 15397

CAS Registry No.: 577784-91-9

Formal Name: 2-chloro-N-[3-(5-chloro-3-methylbenzo[b]thien-2-yl)-1-methyl-1H-pyrazol-5-yl]-benzenesulfonamide

MF: C₁₉H₁₅Cl₂N₃O₂S₂

FW: 452.4

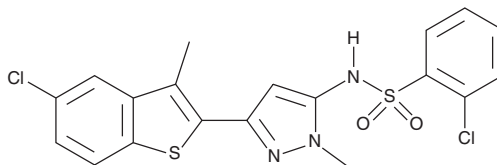
Purity: ≥98%

UV/Vis.: λ_{max}: 243, 299 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: As supplied, 2 years from the QC date provided on the Certificate of Analysis, when stored properly



Laboratory Procedures

ML-60218 is supplied as a crystalline solid. A stock solution may be made by dissolving the ML-60218 in the solvent of choice. ML-60218 is soluble in DMSO at a concentration of approximately 10 mg/ml.

Description

ML-60218 is a small molecule inhibitor of RNA polymerase III in both *S. cerevisiae* and human with IC₅₀ values of 32 and 27 μM, respectively.¹ At 200 μM, it has been shown to inhibit 88, 94, and 90% of the *in vitro* transcription activity of *S. cerevisiae*, *C. albicans*, and human RNA polymerase III, respectively. In combination with HDAC inhibition, ML-60218 at 100 μM has been used to target transcription during exocrine pancreatic epithelial proliferation in a zebrafish model of pancreatic cancer.²

References

1. Wu, L., Pan, J., Thoroddsen, V., *et al.* Novel small-molecule inhibitors of RNA polymerase III. *Eukaryot. Cell* **2**(2), 256-264 (2003).
2. Yee, N.S., Zhou, W., Chun, S.G., *et al.* Targeting developmental regulators of zebrafish exocrine pancreas as a therapeutic approach in human pancreatic cancer. *Biol. Open* **1**(4), 295-307 (2012).

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, 04/27/2016

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