

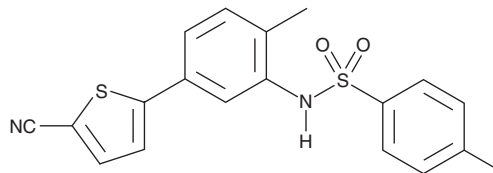
# PRODUCT INFORMATION



**ELR510444**

Item No. 17696

**CAS Registry No.:** 1233948-35-0  
**Formal Name:** N-[5-(5-cyano-2-thienyl)-2-methylphenyl]-4-methyl-benzenesulfonamide  
**MF:** C<sub>19</sub>H<sub>16</sub>N<sub>2</sub>O<sub>2</sub>S<sub>2</sub>  
**FW:** 368.5  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 309 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

ELR510444 is supplied as a crystalline solid. A stock solution may be made by dissolving the ELR510444 in the solvent of choice, which should be purged with an inert gas. ELR510444 is soluble in the organic solvent methyl acetate at a concentration of approximately 10 mg/ml.

## Description

ELR510444 is an inhibitor of tubulin polymerization (IC<sub>50</sub> = 10 μM) that binds to tubulin at the colchicine (Item No. 9000760) binding site.<sup>1</sup> It induces microtubule depolymerization in A-10 cells (EC<sub>50</sub> = 21 nM). ELR510444 inhibits growth in a cancer cell line panel (IC<sub>50</sub>s = 9-43 nM) via formation of multiple spindles and induction of mitotic arrest. Oral administration of ELR510444 (3-6 mg/kg) reduces tumor size in an MDA-MB-231 breast cancer mouse xenograft model in a dose-dependent manner. ELR510444 also inhibits hypoxia-inducible factor 1α (HIF-1α) in RCC4 cells in a concentration-dependent manner.<sup>2</sup>

## References

1. Risinger, A.L., Westbrook, C.D., Encinas, A., *et al.* ELR510444, a novel microtubule disruptor with multiple mechanisms of action. *J. Pharmacol. Exp. Ther.* **336**(3), 652-660 (2011).
2. Carew, J.S., Esquivel, J.A., II, Espitia, C.M., *et al.* ELR510444 inhibits tumor growth and angiogenesis by abrogating HIF activity and disrupting microtubules in renal cell carcinoma. *PLoS One* **7**(1), e31120 (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, 12/14/2022

## 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