

# PRODUCT INFORMATION



## FOX M1 Inhibitor III

Item No. 36932

**CAS Registry No.:** 313380-27-7  
**Formal Name:** 3-amino-N-(4-fluorophenyl)-6-(2-thienyl)-4-(trifluoromethyl)-thieno[2,3-b]pyridine-2-carboxamide

**Synonyms:** FDI-6,  
Forkhead Domain Inhibitory Compound 6,  
NCGC00099374

**MF:** C<sub>19</sub>H<sub>11</sub>F<sub>4</sub>N<sub>3</sub>OS<sub>2</sub>

**FW:** 437.4

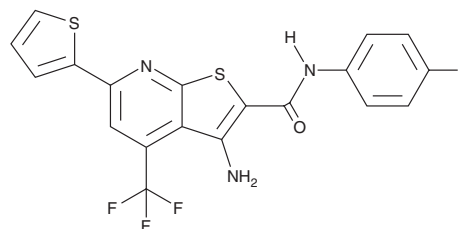
**Purity:** ≥98%

**UV/Vis.:** λ<sub>max</sub>: 285, 327, 354, 369, 410 nm

**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

FOX M1 inhibitor III is supplied as a solid. A stock solution may be made by dissolving the FOX M1 inhibitor III in the solvent of choice, which should be purged with an inert gas. FOX M1 inhibitor III is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of FOX M1 inhibitor III in these solvents is approximately 1 mg/ml.

### Description

FOX M1 inhibitor III is an inhibitor of the DNA-binding activity of forkhead box protein M1 (FOX M1; IC<sub>50</sub> = 22.5 μM).<sup>1</sup> It binds to FOX M1 and is selective for inhibiting FOX M1 transcriptional activity over FOXA1, FOXA2, FXP2, and GATA-binding factor 1 (GATA-1) transcriptional activity. FOX M1 inhibitor III inhibits the growth of MCF-7 and MDA-MB-231 breast and PEO1 ovarian cancer cells (GI<sub>50</sub>s = 18, 21.8, and 18.1 μM, respectively).

### Reference

1. Gormally, M.V., Dexheimer, T.S., Marsico, G., *et al.* Suppression of the FOX M1 transcriptional programme via novel small molecule inhibition. *Nat. Commun.* **5**, 5165 (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, 11/17/2022

#### 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