

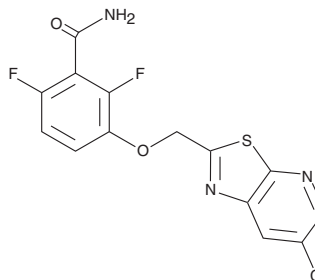
PRODUCT INFORMATION



PC190723

Item No. 13518

CAS Registry No.: 951120-33-5
Formal Name: 3-[[6-chlorothiazolo[5,4-b]pyridin-2-yl)methoxy]-2,6-difluoro-benzamide
MF: C₁₄H₈ClF₂N₃O₂S
FW: 355.0
Purity: ≥98%
UV/Vis.: λ_{max}: 218, 298 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

PC190723 is supplied as a crystalline solid. A stock solution may be made by dissolving the PC190723 in the solvent of choice, which should be purged with an inert gas. PC190723 is soluble (≥10 mg/ml) in DMSO.

Description

PC190723 is an inhibitor of bacterial FtsZ, a homolog of mammalian β-tubulin.¹ It is active against *B. subtilis*, *S. epidermidis*, *S. hemolyticus*, and *S. aureus*, including methicillin-resistant *S. aureus* (MRSA) and multidrug-resistant *S. aureus* (MDRSA; MICs = 0.5-1 μg/ml). PC190723 (10 and 30 mg/kg) increases survival in a mouse model of lethal *S. aureus* infection.¹

Reference

1. Haydon, D.J., Stokes, N.R., Ure, R., *et al.* An inhibitor of FtsZ with potent and selective anti-staphylococcal activity. *Science* **321**, 1673-1675 (2008).

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, 05/02/2025

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