

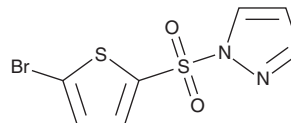
PRODUCT INFORMATION



Qstatin

Item No. 36609

CAS Registry No.: 902688-24-8
Formal Name: 1-[(5-bromo-2-thienyl)sulfonyl]-1H-pyrazole
MF: C₇H₅BrN₂O₂S₂
FW: 293.2
Purity: ≥98%
UV/Vis.: λ_{max}: 230, 252, 275 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

Qstatin is supplied as a solid. A stock solution may be made by dissolving the qstatin in the solvent of choice, which should be purged with an inert gas. Qstatin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of qstatin in these solvents is approximately 1, 5, and 3 mg/ml, respectively.

Description

Qstatin is an inhibitor of quorum sensing.¹ It inhibits activation of the quorum-sensing regulator LuxR in a β-galactosidase reporter assay when used at a concentration of 25 μM.

Reference

1. Dorn, S.K., Newman, J.D., van Kessel, J.C., *et al.* Synthesis and biological assay of small-molecule quorum sensing inhibitors: A three-week course-based undergraduate research experience. *J. Chem. Educ.* **98**(11), 3533–3541 (2021).

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, 10/05/2022

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