

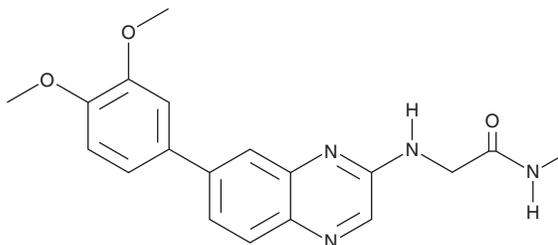
PRODUCT INFORMATION



BQR-695

Item No. 27482

CAS Registry No.: 1513879-21-4
Formal Name: 2-[[7-(3,4-dimethoxyphenyl)-2-quinoxaliny]amino]-N-methylacetamide
MF: C₁₉H₂₀N₄O₃
FW: 352.4
Purity: ≥98%
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

BQR-695 is supplied as a solid. A stock solution may be made by dissolving the BQR-695 in the solvent of choice, which should be purged with an inert gas. BQR-695 is soluble in the organic solvent DMSO.

Description

BQR-695 is an antimalarial compound.¹ It inhibits recombinant *P. vivax* and human phosphatidylinositol 4-kinase (PI4K; IC₅₀s = 3.5 and 90 nM, respectively). BQR-695 induces schizont-stage arrest and death of *P. falciparum* (IC₅₀ = ~70 nM) without inducing cytotoxicity in mature red blood cells (RBCs).

Reference

1. McNamara, C.W., Lee, M.C., Lim, C.S., *et al.* Targeting Plasmodium phosphatidylinositol 4-kinase to eliminate malaria. *Nature* **504(7479)**, 248-253 (2013).

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, 11/08/2022

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