

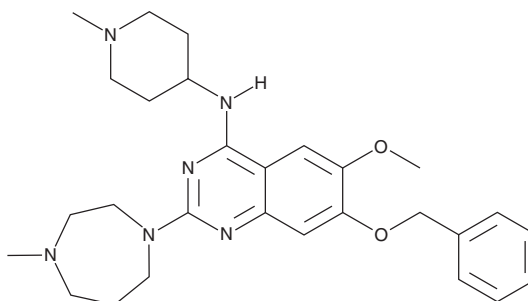
PRODUCT INFORMATION



TM2-115

Item No. 42901

CAS Registry No.: 1197196-47-6
Formal Name: 2-(hexahydro-4-methyl-1H-1,4-diazepin-1-yl)-6-methoxy-N-(1-methyl-4-piperidiny)-7-(phenylmethoxy)-4-quinazolinamine
MF: C₂₈H₃₈N₆O₂
FW: 490.7
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

TM2-115 is supplied as a solid. A stock solution may be made by dissolving the TM2-115 in the solvent of choice, which should be purged with an inert gas. TM2-115 is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

TM2-115 is an inhibitor of *P. falciparum* histone methyltransferase (HMT; IC₅₀ = 43 nM).¹ It is selective for *P. falciparum* HMT over the human HMT G9a (IC₅₀ = >1,000 nM). TM2-115 (1 μM) decreases trimethylation of lysine 4 on histone H3 (H3K4me3) in *P. falciparum*.² It is active against chloroquine-resistant and -sensitive strains of *P. falciparum* (IC₅₀s = 100 and 116 nM, respectively). TM2-115 (1 μM) decreases parasite levels in *P. falciparum*-infected isolated human erythrocytes. It reduces parasitemia in *P. berghei*-infected mice when administered at a dose of 40 mg/kg.

References

1. Sundriyal, S., Chen, P.B., Lubin, A.S., *et al.* Histone lysine methyltransferase structure activity relationships that allow for segregation of G9a inhibition and anti-Plasmodium activity. *Medchemcomm* **8**(5), 1069-1092 (2017).
2. Malmquist, N.A., Moss, T.A., Mecheri, S., *et al.* Small-molecule histone methyltransferase inhibitors display rapid antimalarial activity against all blood stage forms in Plasmodium falciparum. *Proc. Natl. Acad. Sci. USA* **109**(41), 16708-16713 (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, 03/04/2025

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