

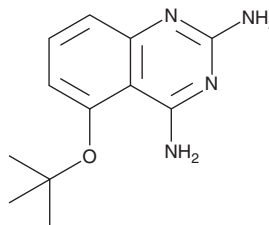
PRODUCT INFORMATION



HZ-1157

Item No. 23436

CAS Registry No.: 1009734-33-1
Formal Name: 5-(1,1-dimethylethoxy)-2,4-quinazolinediamine
MF: $C_{12}H_{16}N_4O$
FW: 232.3
Purity: $\geq 98\%$
UV/Vis.: λ_{\max} : 240, 335 nm
Supplied as: A crystalline 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

HZ-1157 is supplied as a crystalline solid. A stock solution may be made by dissolving the HZ-1157 in the solvent of choice, which should be purged with an inert gas. HZ-1157 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of HZ-1157 in ethanol is approximately 1 mg/ml and approximately 5 mg/ml in DMSO and DMF.

Description

HZ-1157 is an inhibitor of hepatitis C virus (HCV) nonstructural protein 3/4A (NS3/4A) protease ($IC_{50} = 1 \mu\text{M}$ in an NS3/4A secreted embryonic alkaline phosphatase assay).¹ It inhibits infectious HCV virus replication in Huh7.51 cells ($IC_{50} = 0.82 \mu\text{M}$). It also inhibits dengue virus serotype 2 replicon activity ($EC_{50} = 2.8 \text{ nM}$ in BHK-D2RepT cells with high NS3 levels) without exhibiting cytotoxicity ($CC_{50} = >10 \mu\text{M}$).²

References

1. Yu, Y., Jing, J.F., Tong, X.K., *et al.* Discovering novel anti-HCV compounds with inhibitory activities toward HCV NS3/4A protease. *Acta. Pharmacol. Sin.* **35(8)**, 1074-1081 (2014).
2. Chao, B., Tong, X.K., Tang, W., *et al.* Discovery and optimization of 2,4-diaminoquinazoline derivatives as a new class of potent dengue virus inhibitors. *J. Med. Chem.* **55(7)**, 3135-3143 (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, 11/15/2022

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