

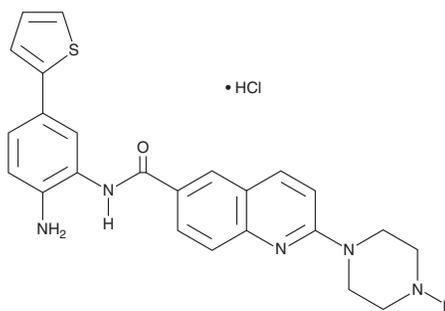
# PRODUCT INFORMATION



## ACY-957 (hydrochloride)

Item No. 43477

**Formal Name:** N-[2-amino-5-(2-thienyl)phenyl]-2-(1-piperazinyl)-6-quinolinecarboxamide, monohydrochloride  
**MF:** C<sub>24</sub>H<sub>23</sub>N<sub>5</sub>OS • HCl  
**FW:** 466.0  
**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

ACY-957 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the ACY-957 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. ACY-957 (hydrochloride) is sparingly soluble (1-10 mg/ml) in DMSO.

### Description

ACY-957 is an inhibitor of histone deacetylase 1 (HDAC1) and HDAC2 (IC<sub>50</sub>s = 7 and 18 nM, respectively).<sup>1</sup> It is selective for HDAC1 and HDAC2 over HDAC3 (IC<sub>50</sub> = 1,300 nM) and does not inhibit HDAC4-9 at concentrations up to 20 μM. ACY-957 (1 μM) induces expression of *HBG*, the gene encoding γ-globin, in peripheral blood mononuclear cells (PBMCs) isolated from patients with sickle cell disease.

### Reference

1. Shearstone, J.R., Golonzhka, O., Chonkar, A., *et al.* Chemical inhibition of histone deacetylases 1 and 2 induces fetal hemoglobin through activation of GATA2. *PLoS One* **11**(4), e0153767 (2016).

#### 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, 07/31/2025

#### 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