

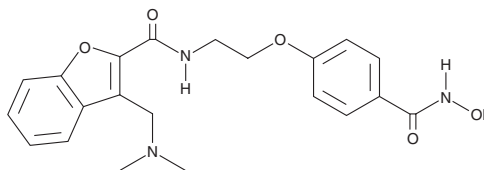
# PRODUCT INFORMATION



## PCI 24781

Item No. 20059

**CAS Registry No.:** 783355-60-2  
**Formal Name:** 3-[[dimethylamino)methyl]-N-[2-[4-[(hydroxyamino)carbonyl]phenoxy]ethyl]-2-benzofurancarboxamide  
**Synonyms:** CRA-024781, CRA-24781  
**MF:** C<sub>21</sub>H<sub>23</sub>N<sub>3</sub>O<sub>5</sub>  
**FW:** 397.4  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 265 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

PCI 24781 is supplied as a crystalline solid. A stock solution may be made by dissolving the PCI 24781 in the solvent of choice, which should be purged with an inert gas. PCI 24781 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of PCI 24781 in these solvents is approximately 3, 30, and 50 mg/ml, respectively.

PCI 24781 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, PCI 24781 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. PCI 24781 has a solubility of approximately 0.1 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

PCI 24781 is an HDAC inhibitor that predominantly targets HDAC1 with a K<sub>i</sub> value of 7 nM.<sup>1</sup> It inhibits HDAC2, HDAC3/SMRT, HDAC6, HDAC8, and HDAC10 with K<sub>i</sub> values of 19, 8.2, 17, 280, and 24 nM, respectively.<sup>1</sup> PCI 24781 can prevent the growth of tumor cells *in vitro*, inducing apoptosis, as well as in mice harboring HCT116 or DLD-1 colon tumor xenografts.<sup>1</sup>

### Reference

1. Buggy, J.J., Cao, Z.A., Bass, K.E., *et al.* CRA-024781: A novel synthetic inhibitor of histone deacetylase enzymes with antitumor activity *in vitro* and *in vivo*. *Mol. Cancer Ther.* **5**(5), 1309-1317 (2006).

#### 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, 12/15/2022

#### 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