

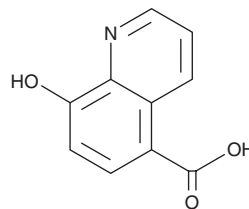
# PRODUCT INFORMATION



## IOX1

Item No. 11572

CAS Registry No.: 5852-78-8  
Formal Name: 8-hydroxy-5-quinolinecarboxylic acid  
MF:  $C_{10}H_7NO_3$   
FW: 189.2  
Purity:  $\geq 97\%$   
UV/Vis.:  $\lambda_{max}$ : 241, 320 nm  
Supplied as: A crystalline solid  
Storage:  $-20^{\circ}C$   
Stability:  $\geq 4$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

IOX1 is supplied as a crystalline solid. A stock solution may be made by dissolving the IOX1 in the solvent of choice, which should be purged with an inert gas. IOX1 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of IOX1 in these solvents is approximately 20 and 10 mg/ml, respectively.

IOX1 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, IOX1 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. IOX1 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

IOX1 is a broad-spectrum inhibitor of 2OG oxygenases that does not require application in a pro-drug formulation. IOX1 inhibits JMJD2A, JMJD2E and the 2OG oxygenases PHF8, PHD2, and FIH with  $IC_{50}$  values of 1.7, 2.4, 13.3, 14.3, and 20.5  $\mu M$ , respectively.<sup>1</sup> IOX1 inhibits H3K9me3 demethylation by JMJD2A in HeLa cells with an  $IC_{50}$  value of 87  $\mu M$ .<sup>1</sup>

### Reference

1. King, O.N.F., Li, X.S., Sakurai, M., *et al.* Quantitative high-throughput screening identifies 8-hydroxyquinolines as cell-active histone demethylase inhibitors. *PLoS One* **5(11)**, 1-12 (2010).

#### 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, 11/10/2022

#### 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