

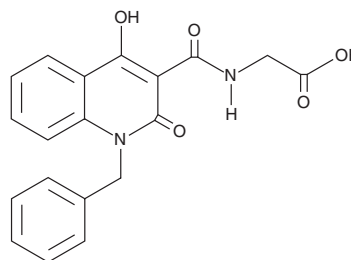
# PRODUCT INFORMATION



## IOX2

Item No. 11573

**CAS Registry No.:** 931398-72-0  
**Formal Name:** N-[[1,2-dihydro-4-hydroxy-2-oxo-1-(phenylmethyl)-3-quinoliny]carbonyl]-glycine  
**MF:** C<sub>19</sub>H<sub>16</sub>N<sub>2</sub>O<sub>5</sub>  
**FW:** 352.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 232, 289, 331 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

IOX2 is supplied as a crystalline solid. A stock solution may be made by dissolving the IOX2 in the solvent of choice, which should be purged with an inert gas. IOX2 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of IOX2 in these solvents is approximately 5 and 14 mg/ml, respectively.

IOX2 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, IOX2 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. IOX2 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

IOX2 is an inhibitor of hypoxia-inducible factor prolyl hydroxylase 2 (HIF-PH2; IC<sub>50</sub> = 21 nM) with over 100-fold selectivity compared to inhibition of JMJD2A, JMJD2C, JMJD2E, JMJD3, or the 2OG oxygenase FIH (IC<sub>50</sub>s > 100 μM).<sup>1</sup> IOX2 is active in cells, inhibiting HIF-1α hydroxylation in RCC4 cells at 50 μM.<sup>1</sup> See the Structural Genomics Consortium (SGC) website for more information.

### Reference

1. IOX2: A selective inhibitor of the hypoxia inducible factor (HIF) prolyl-hydroxylases. Retrieved June 13, 2012, from [http://www.thesgc.org/scientists/chemical\\_probes/IOX2](http://www.thesgc.org/scientists/chemical_probes/IOX2).

#### 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, 11/17/2022

#### 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](http://WWW.CAYMANCHEM.COM)