

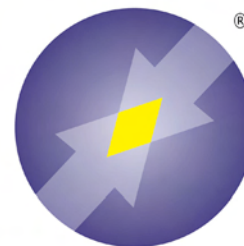
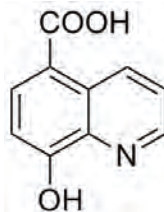
# IOX1

Catalog Number P022-5MG

Catalog Number P022-25MG

## FEATURES

- Inhibitor of *Jumonji* HDMs
- Broad spectrum 2-oxoglutarate oxygenase inhibitor
- Cell Permeable



ARBOR  
ASSAYS

## INTRODUCTION

Inhibits the JMJD family of 2-oxoglutarate-dependent histone demethylases. IC<sub>50</sub> = 0.12, 0.07, 0.2, 0.3, 0.6, and 1 μM for JMJD3, JMJD1A, JMJD2A, JMJD2E, JMJD2C and UTX respectively<sup>1</sup>. Is a broad spectrum 2-oxoglutarate oxygenase inhibitor which may be used to study hypoxic signaling<sup>1</sup>. Cell permeable.

<b>FORM:</b>	Yellow powder
<b>MOLECULAR WEIGHT:</b>	189.2
<b>STORAGE:</b>	Room temperature desiccated
<b>FORMULA:</b>	C <sub>10</sub> H <sub>7</sub> NO <sub>3</sub>
<b>CAS NUMBER:</b>	5852-78-8
<b>OTHER NAMES:</b>	8-Hydroxy-5-quinolinecarboxylic acid; (5-Carboxy-8HQ)
<b>USES:</b>	Soluble to 20 mg/mL in DMSO

## REFERENCES:

1. ON King et al. PLoS One 2010 5: e15535

**FOR RESEARCH USE ONLY**