

Catalog No: 14056, 14057

Format: 25 mg, 5 mg

Chemical Properties:

MW = 189.2 $C_{10}H_7NO_3$

CAS 5852-78-8

Physical Properties: Yellow powder

Names: IOX-1, 8-Hydroxy-5-quinolinecarboxylic acid; 5-Carboxy-8HQ

Pharmacology: Inhibits the JMJD family of 2-oxoglutarate-dependent histone demethylases. IC_{50} = 0.12, 0.07, 0.2, 0.3, 0.6 and 1 M for JMJD3, JMJD1A, JMJD2A, JMJD2E, JMJD2C and UTX respectively (ref 1). A broad spectrum 2-oxoglutarate oxygenase inhibitor which may be used to study hypoxic signaling (ref 1). Cell permeable.

Solubilization: May be dissolved in DMSO (20 mg/ml)

Fluorescent Properties: N/A

Quality Control:

>98% (HPLC); NMR (Conforms)

References:

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

Storage and Guarantee: Store desiccated as supplied at room temperature for up to 1 year. Store solutions at -20°C for up to 3 months. This product is guaranteed for 6 months from date of arrival.

