

NSC-663284

Catalog No: 14122, 14123

Format: 25 mg, 5 mg

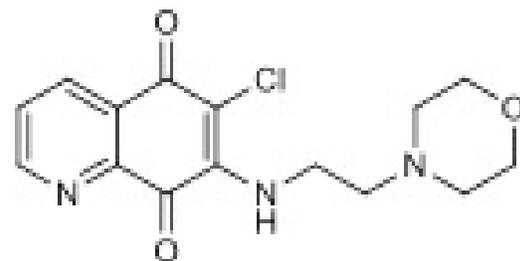
**Chemical Properties:**

MW = 321.8

C<sub>15</sub>H<sub>16</sub>ClN<sub>3</sub>O<sub>3</sub>

CAS 383907-43-5

Physical Properties: Red powder

**Names:** NSC-663284, 6-Chloro-7-[[2-(4-morpholinyl)ethyl]amino]-5,8-quinolinedione

**Pharmacology:** Potent and selective inhibitor of the protein lysine methyltransferase SETD8 (IC<sub>50</sub> = 0.21 μM) which efficiently suppresses cellular H4K20me1 at doses lower than 5 μM (ref 1). Also inhibits CDC25 phosphatases (ref 2). Arrests cells at G<sub>1</sub> and G<sub>2</sub>/M and inhibits cdk1 and 2 activation (ref 3). Inhibits the growth of a variety of human tumor cell lines (ref 4).

**Solubilization:** May be dissolved in DMSO (15 mg/ml); or ethanol (55 mg/ml)

**Fluorescent Properties:** N/A

**Quality Control:**

>98% (TLC); NMR (Conforms)

## References:

1. G Blum *et al.* *ACS Chem. Biol.* 2014, 9:2471
2. JS Lazo *et al.* *J. Med. Chem.* 2001, 44:4042
3. L Pu *et al.* *J. Biol. Chem.* 2002, 277:46877
4. Y Han *et al.* *J. Pharmacol. Exp. Ther.* 2004, 309:64

**Storage and Guarantee:** Store desiccated as supplied at -20° for up to 2 years. Store solutions at -20°C for up to 1 month. This product is for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.