

Catalog No: 14104, 14105

Format: 25 mg, 5 mg

Chemical Properties:

MW = 334.3 $C_{19}H_{14}N_2O_4$

CAS 48208-26-0

Physical Properties: Tan powder, MP = 104°C

Names: RG-108, N-Phthalyl-L-tryptophan; or (S)-2-(1,3-dioxoisoindolin-2-yl)

-3-(1H-indol-3-yl)propanoic acid

Pharmacology: A non-nucleoside DNA methyltransferase inhibitor (IC_{50} = 115 nM)(ref 1). Inhibits DNA methylation in human and tumor cell lines without toxicity (ref 2). Enhances the reversion of neural progenitor cells to the pluripotent state (ref 3).

Solubilization: May be dissolved in DMSO (100 mg/ml); Ethanol (100 mg/ml)

Fluorescent Properties: N/A

Quality Control:

>98% (TLC); NMR (Conforms)

References:

1. B Brueckner et al. Cancer Res. 2005, 65:6305

2. C Stresemann et al. Cancer Res. 2006, 66:2794

3. I Szablowska-Gadomska et al. Acta. Neurobiol. Exp. 2012, 72:337

Storage and Guarantee: Store desiccated as supplied at room temperature for up to 2 years. Store solutions at -20°C for up to 1 month. This product is guaranteed for 6 months from date of arrival.