

RG-108

Epigenetic Enzyme Inhibitor

Catalog # D353-904C

Lot # P3320-31

CAS # 48208-26-0

Product Description

Molecular Formula: C₁₉H₁₄N₂O₄

Appearance: Tan powder

Melting Point: 104°C

Molecular Weight: 334.3

Purity: >98% (TLC); NMR (Conforms)

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

Alias

N-Phthalyl-L-tryptophan

Specific Activity

DNA Methyltransferase Inhibitor

Storage and Stability

Store desiccated as supplied at ambient temperature for up to 2 years. Store solutions at -20°C for up to 1 month.

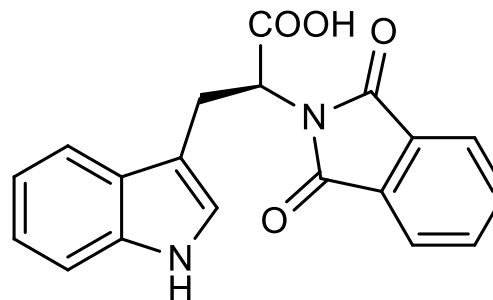
Scientific Background

A non-nucleoside DNA methyltransferase inhibitor. IC₅₀=115 nM¹ Inhibits DNA methylation in human tumor cell lines without toxicity². Enhances the reversion of neural progenitor cells to the pluripotent state³.

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

Molecular Structure



RG-108

Epigenetic Enzyme Inhibitor

Catalog # D353-904C

CAS # 48208-26-0

Lot # P3320-31

Purity >98% (TLC)

Format Tan powder

Solubilization May be dissolved in DMSO (100 mg/ml); or ethanol (100 mg/ml)

Storage & Shipping Store desiccated as supplied at ambient temperature for up to 2 years. Store solutions at -20°C for up to 1 month.