
RG 108

(Catalog No. M25000)

Alternative Names

N-Phthalyl-L-tryptophan

Description

Non-nucleoside DNA methyltransferase inhibitor that blocks the enzyme active site. In vitro inhibits DNA methylation in human cancer cell lines without detectable toxicity. Demethylates and reactivates epigenetically silenced tumor suppressor genes.

CAS Number

48208-26-0

Chemical Formula

$C_{19}H_{14}N_2O_4$

Molecular Weight

334.33

Size

10 mg

Purity

≥98%

Solubility

Soluble to 100 mM in DMSO and to 100 mM in Ethanol

Storage Temp

-20°C

Ordering Information:

Products	Size	Cat. No.
RG 108	10 mg	M25000 -1

This product is for research purposes only. Not intended for use in diagnostic procedures.