

Catalog No: 14048, 14049

Format: 25 mg, 5 mg

Chemical Properties:

MW = 339.4 $C_{20}H_{25}N_3O_2$

CAS 537034-17-6

Physical Properties: Tan powder, MP = 168-170°C

Names: BML 210, N1-(2-aminophenyl)-N8-phenyloctanediamide

Pharmacology: A novel non-hydroxamic HDAC inhibitor. ($IC_{50} = 5-10 \text{ M}$) (HeLa nuclear extract). Induces growth inhibition, proapoptotic and differentiation effects on human leukemia cell lines.

Solubilization: May be dissolved in DMSO (25 mg/ml); Ethanol (10 mg/ml,

warm)

Fluorescent Properties: N/A

Quality Control:

98% (TLC); NMR (Conforms)

References:

1. J Savickiene et al. Eur. J. Pharmacol. 2006, 549:9

2. D Herman et al. Nature Chem. Biol. 2006, 10:551

Storage and Guarantee: Store desiccated as supplied at 4°C for up to 2 years. Store solutions at -20°C for up to 3 months. This product is guaranteed for 6 months from date of arrival.