

BML-210

Epigenetic Enzyme Inhibitor

Catalog # H83-904C

Lot # P3320-6

CAS # 537034-17-6

Product Description

Molecular Formula: C₂₀H₂₅N₃O₂

Appearance: Tan powder

Melting Point: 168-170°C

Molecular Weight: 339.4

Purity: 98% ; NMR (Conforms)

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

Alias

N1-(2-aminophenyl)-N8-phenyloctanediamide

Specific Activity

HDAC Inhibitor

Storage and Stability

Store desiccated as supplied at 4°C for up to 2 years.

Store solutions at -20°C for up to 3 months

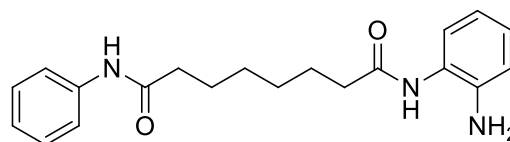
Scientific Background

A novel non-hydroxamic acid HDAC inhibitor. IC₅₀=5-10 μM (HeLa nuclear extract). Induces growth inhibition, proapoptotic and differentiation effects on human leukemia cell lines.

References

1. J. Savickiene et al. Eur. J. Pharmacol. 2006 549 9
2. D. Herman et al. Nature Chem. Biol. 2006 10 551

Molecular Structure



BML-210

Epigenetic Enzyme Inhibitor

| | |
|--------------------|--|
| Catalog # | H83-904C |
| CAS # | 537034-17-6 |
| Lot # | P3320-6 |
| Purity | 98% |
| Format | Tan powder |
| Solubilization | May be dissolved in DMSO (25 mg/ml); or ethanol (10 mg/ml, warm) |
| Storage & Shipping | Store desiccated as supplied at 4°C for up to 2 years. Store solutions at -20°C for up to 3 months |

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.