## BML-278



Catalog No: 14024, 14025 Format: 25 mg, 5 mg

## **Chemical Properties:**

 $\label{eq:mw} \begin{array}{l} \mathsf{MW} = 391.5\\ \mathsf{C}_{24}\mathsf{H}_{25}\mathsf{NO}_4\\ \\ \mathsf{CAS} \text{ Not found}\\ \\ \mathsf{Physical Properties: off white powder, MP = 141-143^\circ\mathsf{C}} \end{array}$ 

Names: BML-278, N-Benzyl-3,5-dicarbethoxy-4-phenyl-1,4-dihydropyridine

**Pharmacology:** A structurally novel activator of SIRT1 ( $EC_{150}$ = 1 M). Displays selectivity over SIRT2 and SIRT3 ( $EC_{150}$ = 25 M and 50 M respectively). Cell permeable.

Solubilization: May be dissolved in DMSO (5 mg/ml)

Fluorescent Properties: N/A

Quality Control: >98% (TLC); NMR (Conforms)

References: 1. A Mai *et al. J. Med. Chem.* 2009, 52:5496

**Storage and Guarantee:** Store desiccated as supplied at room temperature 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.

