

## BML-278

Epigenetic Enzyme Activator

**Catalog # S35-904C**

Lot # P3320-7

CAS # N/F

### Product Description

Molecular Formula: C<sub>24</sub>H<sub>25</sub>NO<sub>4</sub>

Appearance: Off white powder

Melting Point: 141-143°C

Molecular Weight: 391.5

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

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

### Alias

N-Benzyl-3,5-dicarbethoxy-4-phenyl-1,4-dihydropyridine

### Specific Activity

SIRT1 Activator

### Storage and Stability

Store desiccated as supplied at room temperature for up to 2 years. Store solutions at -20°C for up to 3 months.

### Scientific Background

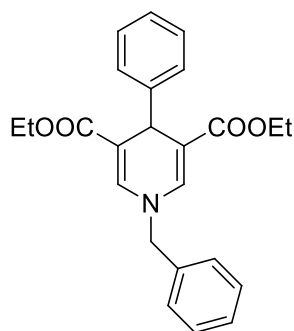
A structurally novel activator of SIRT1 (EC<sub>150</sub>=1 μM).

Displays selectivity over SIRT2 and 3 (EC<sub>150</sub>=25 and 50 μM respectively). Cell permeable.

### References

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

### Molecular Structure



## BML-278

Epigenetic Enzyme Activator

Catalog # S35-904C

CAS # N/F

Lot # P3320-7

Purity >98% (TLC)

Format Off white powder

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

Storage & Shipping Store desiccated as supplied at room temperature 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](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

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