

BML-278

SIRT activator

Novel Sirt activator structurally based on a dihydropyridine scaffold. EC_{150} (effective concentration able to increase enzyme activity by 150%) = $1\mu M$ for Sirt1 with weaker effect on Sirt2 and Sirt3 (25 and $50\mu M$ respectively). The compound is active in whole cells and arrests cell cycle at G1/S phase in U937 cells.

Citations: 1

[View Online »](#)

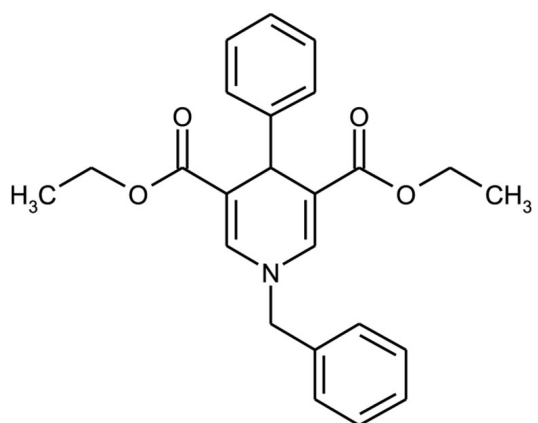
Ordering Information

[Order Online »](#)

BML-GR359-0005	5mg
BML-GR359-0025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at room temperature. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	N-Benzyl-3,5-dicarbethoxy-4-phenyl-1,4-dihydropyridine
Appearance	Off-white solid.
CAS	15301-69-6
Couple Target	SIRT
Couple Type	Activator
Formula	$C_{24}H_{25}NO_4$
MW	391.5
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (5mg/ml).
Source	Synthetic.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)