

BML-266

SIRT inhibitor

Structurally novel SIRT2 inhibitor (IC_{50} =56.7 μ M).

Citations: 1

[View Online »](#)

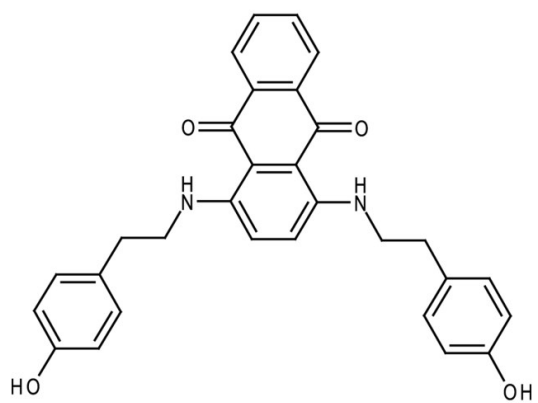
Ordering Information

[Order Online »](#)

BML-GR346-0010	10mg
BML-GR346-0050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 1,4-Di[(4-hydroxyphenethyl)amino]-9,10-dihydroanthracene-9,10-dione

Appearance Dark purple solid.

CAS 96969-83-4

Couple Target SIRT, SIRT2

Couple Type Inhibitor

Formula $C_{30}H_{26}N_2O_4$

Identity Determined by NMR.

MW 478.5

Purity ≥95% (TLC)

Solubility Soluble in DMSO (25mg/ml).

Last modified: May 29, 2024



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-
fr@enzolifesciences.com](mailto:info-fr@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)