

Mirin

MRN inhibitor

The MRN-ATM pathway is essential for sensing the signaling from DNA double-strand breaks. Mirin is an inhibitor of MRN discovered using a forward chemical genetic screen. It prevents MRN-dependent activation of ATM without affecting ATM protein kinase activity and it inhibits Mre11-associated exonuclease activity. $IC_{50}=66\mu M$. It abolishes the G2/M checkpoint and homology-dependent repair in mammalian cells.

Citations: 3

[View Online »](#)

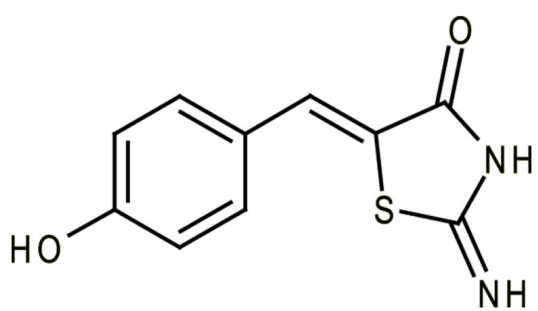
Ordering Information

[Order Online »](#)

BML-GR357-0005	5mg
BML-GR357-0025	25mg

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 Z-5-(4-Hydroxybenzylidene)-2-imino-1,3-thiazolidin-4-one

Appearance Yellow solid.

CAS 299953-00-7

Couple Target MRN

Couple Type Inhibitor

Formula $C_{10}H_8N_2O_2S$

MW 220.3

Purity ≥95% (TLC)

Solubility Soluble in DMSO (25mg/ml) or DMF (25mg/ml).

Source Synthetic.

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)