

1,5-Isoquinolinediol

PARP inhibitor

Potent inhibitor of inducible nitric oxide synthase (iNOS; NOS II) in mouse macrophages. Potent and selective PARP inhibitor. Inhibits nitric oxide (SIN-1) induced PARP activation in rat hepatocytes (at 20µM). Neuroprotectant.

Citations: 10

[View Online »](#)

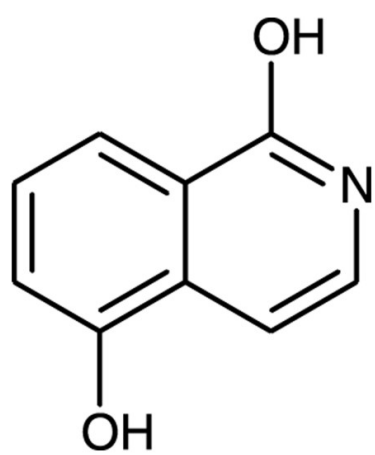
Ordering Information

[Order Online »](#)

BML-AP102-0020	20mg
BML-AP102-0100	100mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Yellow solid.
CAS	5154-02-9
Couple Target	PARP
Couple Type	Inhibitor
Formula	$C_9H_7NO_2$
MW	161.2
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (25mg/ml), dimethyl formamide (25mg/ml) or 100% ethanol (25mg/ml warm). Also soluble in 0.1N NaOH or in methanol; insoluble in water and 0.1N HCl.

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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)