

Dimethyl-bisphenol A

HIF-1 α inhibitor

A HIF-1 α degradation promotor.

Citations: 2

[View Online »](#)

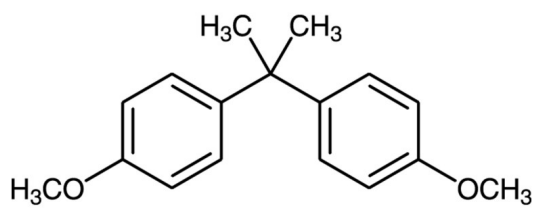
Ordering Information

[Order Online »](#)

BML-GR339-0010	10mg
----------------	------

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 Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance White solid.

CAS 1568-83-8

Couple Target HIF

Couple Type Inhibitor

Formula $C_{17}H_{20}O_2$

MW 256.3

Purity ≥98% (TLC)

Solubility Soluble in DMSO (>25 mg/ml) or ethanol (>25 mg/ml)

Source Synthetic.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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