

Dimethyloxaloylglycine

Prolyl-4-hydroxylase inhibitor

DMOG is a cell permeable prolyl-4-hydroxylase inhibitor which upregulates HIF activity. HIF activation stimulates angiogenesis in several different models. DMOG also inhibits FIH (Factor Inhibiting HIF), an asparaginyl hydroxylase, which enhances the HIF response. It is active *in vivo* and attenuates myocardial injury in a rabbit ischemia reperfusion model (20mg/kg). Is expected to act pro-angiogenic.

Citations: 16

[View Online »](#)

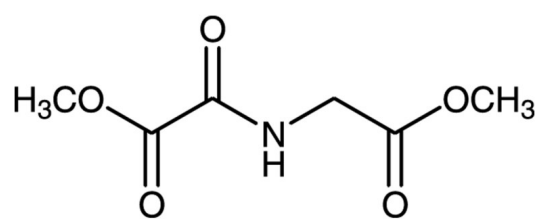
Ordering Information

[Order Online »](#)

BML-EI347-0010	10mg
BML-EI347-0050	50mg

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. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	-20°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	DMOG, N-(Methoxyoxoacetyl)-glycine methyl ester
Appearance	Off-white solid.
CAS	89464-63-1
Couple Target	FIH, HIF, Prolyl-4-hydroxylase
Couple Type	Activator, Inhibitor
Formula	$C_6H_9NO_5$
MW	175.1
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (>25mg/ml), 100% ethanol (>25mg/ml) or dimethyl formamide; soluble also in PBS, pH7.2 (10mg/ml).

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-fr@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