

# 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 »](#)

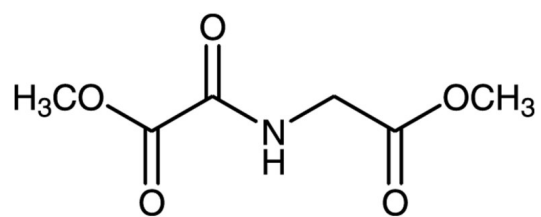
## Ordering Information

[Order Online »](#)

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

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	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.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Dry Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

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



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