

# BML-288

## PDE2 inhibitor

Potent ( $IC_{50}$ =40nM) and selective (290X selectivity over other PDE isoforms) inhibitor of phosphodiesterase Type II (PDE2). Displayed no significant level of binding to neither a collection of 54 receptors and ion channels, nor to a panel of 30 kinases. BML-288 demonstrated no inhibition of 5-lipoxygenase (5-LO) or cyclooxygenase (COX-1). BML-288 is more potent and selective than EHNA (BML-PD129).

Citations: 1

[View Online »](#)

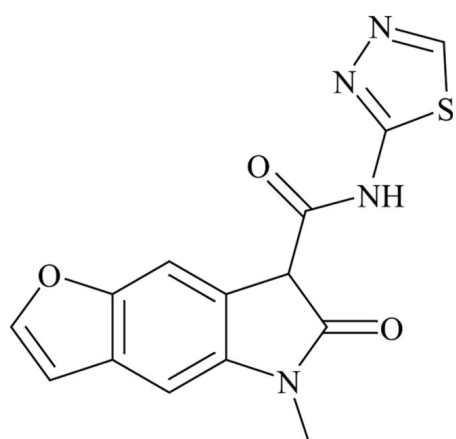
## Ordering Information

[Order Online »](#)

BML-PD200-0025	25mg
----------------	------

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Store under inert gas. Compound will darken on exposure to air. Stable for at least 1 year after the date of receipt when stored at -20°C. Solutions can be stored at -20°C for up to 3 months.
<b>Handling</b>	Solutions can be stored at -20°C for up to 3 months.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	5-Methyl-6-oxo-N-(1,3,4-thiadiazol-2-yl)-7H-furo[2,3-f]indole-7-carboxamide
<b>Appearance</b>	Off-white solid.
<b>Couple Target</b>	PDE
<b>Couple Type</b>	Inhibitor
<b>Formula</b>	$C_{14}H_{10}N_4O_3S$
<b>MW</b>	314.3
<b>MeltingPoint</b>	243-245°C
<b>Purity</b>	≥98% (TLC)
<b>Solubility</b>	Soluble in DMSO (25 mg/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)