

# BRL-50481

PDE inhibitor

The first selective PDE7 inhibitor ( $K_i=180$  nM).

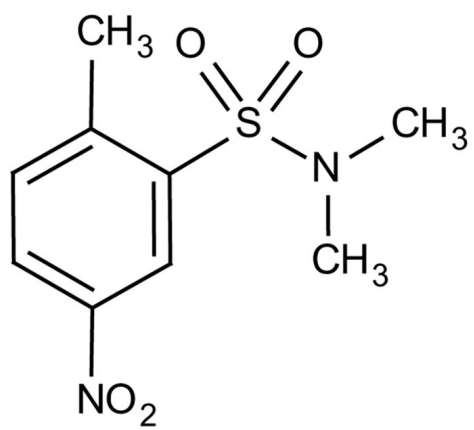
## Ordering Information

[Order Online »](#)

BML-PD120-0050	50mg
----------------	------

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

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

**Regulatory Status** RUO - Research Use Only

## Product Details

Appearance	White to off-white solid.
CAS	433695-36-4
Couple Target	PDE
Couple Type	Inhibitor
Formula	$C_9H_{12}N_2O_4S$
Formulation	Lyophilized.
MW	244.27
MeltingPoint	65-67°C
Purity	≥98% (HPLC)
Source	Synthetic.

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-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)