

Siguazodan

PDE inhibitor

A selective phosphodiesterase PDE III inhibitor, $IC_{50}=0.7\mu M$. Siguazodan is a potent inotropic, vasodilator and antiplatelet agent with sustained duration of action. It synergizes with PDE IV inhibitors (roliprams) in blocking antigen-induced bronchospasm in the guinea pig.

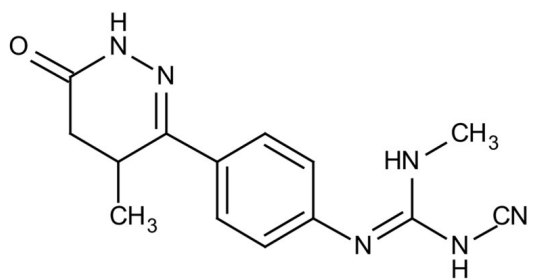
Ordering Information

[Order Online »](#)

BML-PD179-0005	5mg
----------------	-----

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 solid.
CAS	115344-47-3
Couple Target	PDE
Couple Type	Inhibitor
Formula	$C_{14}H_{16}N_6O$
Identity	Determined by 1H-NMR.
MW	284.3
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO.

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)