

Zardaverine

PDE inhibitor

Selective, dual inhibitor of PDE III and PDE IV (IC_{50} s=2.5 and 1.1 μ M respectively). Anti-allergic activity, combined inhibition of PDE III and IV have an additive synergistic effect on inhibition of guinea pig bronchospasm.

Citations: 1

[View Online »](#)

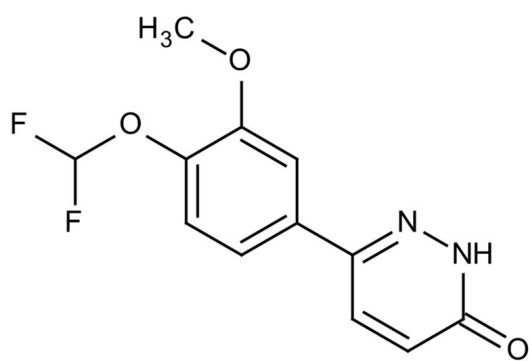
Ordering Information

[Order Online »](#)

BML-PD195-0005	5mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store solutions at -20°C for up to 3 months.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Off-white to yellow solid.
CAS	101975-10-4
Couple Target	PDE
Couple Type	Inhibitor
Formula	$C_{12}H_{10}F_2N_2O_3$
Identity	Determined by 1H -NMR and MS.
MW	268.2
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (25mg/ml).



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