

PGP-4008

P-glycoprotein inhibitor

Selective P-glycoprotein (Pgp) inhibitor. Effectively reverses Pgp mediated drug resistance in NCI/ADR cells, but has no effect on the drug sensitivities of MCF-7/VP cells that overexpress MRP1.

Citations: 2

[View Online »](#)

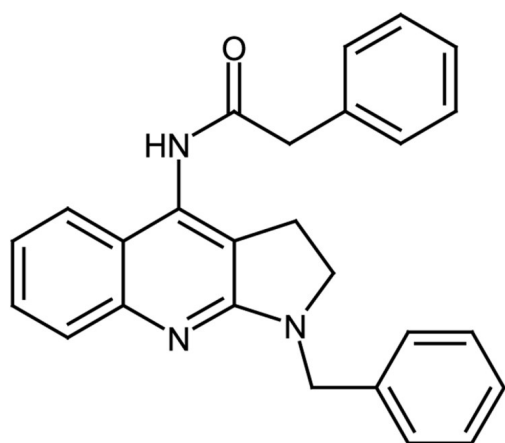
Ordering Information

[Order Online »](#)

ALX-270-290-M002	2mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	N-(1-Benzyl-2,3-dihydro-1H-pyrrolo[2,3-b]quinolin-4-yl)-2-phenylacetamide
Appearance	White to off-white powder.
CAS	365565-02-2
Couple Target	P-glycoprotein
Couple Type	Inhibitor
Formula	$C_{26}H_{23}N_3O$
MW	393.5
Purity	≥93%
Solubility	Soluble in methanol, 100% ethanol (3mg/ml) or DMSO (20mg/ml).

Last modified: May 29, 2024



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-be@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