

ODQ

Guanylyl cyclase inhibitor

ODQ is a potent and selective, direct-acting inhibitor of guanylate cyclase ($IC_{50}=20$ nM). In cerebellar slices ODQ reversibly inhibited cGMP generation in response to endogenous NO or exogenously added NO-donors. ODQ does not inhibit NO-mediated macrophage toxicity (a phenomenon unrelated to cGMP) and does not inhibit particulate guanylyl cyclase or adenylyl cyclase. ODQ is an unique and selective tool to elucidate the physiological importance of the NO-cGMP pathway.

Citations: 19

[View Online »](#)

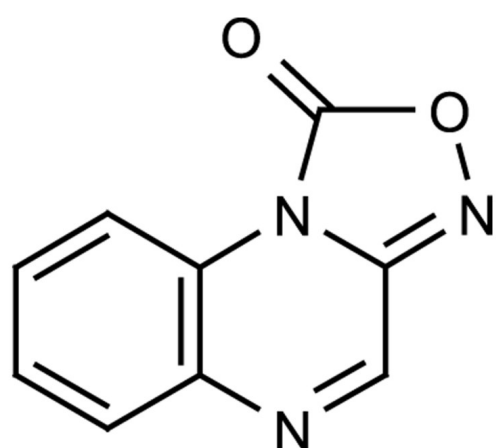
Ordering Information

[Order Online »](#)

ALX-270-034-M050	50mg
------------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Solutions may decompose slowly.
Handling	Protect from light.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Off-white to yellow powder.
CAS	41443-28-1
Couple Target	Guanylyl cyclase
Couple Type	Inhibitor
Formula	$C_9H_5N_3O_2$
MW	187.2
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (5mg/ml), ethyl acetate or 100% ethanol (1.2mg/ml); insoluble in water.

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-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