

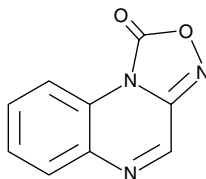
Product Information



ODQ

Item No. 81410

CAS Registry No.: 41443-28-1
Formal Name: 1H-[1,2,4]oxadiazolo[4,3-a]quinoxalin-1-one
MF: C₉H₅N₃O₂
FW: 187.2
Purity: ≥98%
Stability: ≥1 year at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 232, 295 nm



Laboratory Procedures

For long term storage, we suggest that ODQ be stored as supplied at -20°C. It should be stable for at least one year.

ODQ is supplied as a crystalline solid. A stock solution may be made by dissolving the ODQ in an organic solvent purged with an inert gas. ODQ is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of ODQ in these solvents is approximately 100 μM, 15 mg/ml, and 30 mg/ml, respectively.

ODQ is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ODQ should first be dissolved in DMF and then diluted with the aqueous buffer of choice. ODQ has a solubility of approximately 100 μg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

ODQ is a highly selective, irreversible, heme-site inhibitor of soluble guanylyl cyclase.^{1,2} The binding of ODQ is competitive with NO. The inhibition of soluble guanylyl cyclase is time dependent with complete inactivation in 10 minutes at 0.3 μM ODQ.²

References

1. Garthwaite, J., Southam, E., Boulton, C.L. *et al.* Potent and selective inhibition of nitric oxide-sensitive guanylyl cyclase by 1H-[1,2,4]oxadiazolo[4,3-a]quinoxalin-1-one. *Mol. Pharmacol.* **48**, 184-188 (1995).
2. Schrammel, A., Behrends, S., Schmidt, K., *et al.* Characterization of 1H-[1,2,4]oxadiazolo[4,3-a]quinoxalin-1-one as a heme-site inhibitor of nitric oxide-sensitive guanylyl cyclase. *Mol. Pharmacol.* **50**, 1-5 (1996).

Related Products

LY83583 - Item No. 70230 • YC-1 - Item No. 81560 • NS-2028 - Item No. 81600 • Guanylate Cyclase α subunit (soluble) Polyclonal Antibody - Item No. 160895 • Guanylate Cyclase β1 subunit (soluble) Polyclonal Antibody - Item No. 160897

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY; NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants only to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any obligation or liability, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's exclusive remedy and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 05/17/2011