

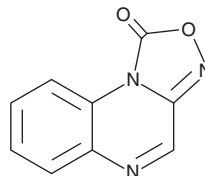
PRODUCT INFORMATION



ODQ

Item No. 81410

CAS Registry No.: 41443-28-1
Formal Name: 1H-[1,2,4]oxadiazolo[4,3-a]quinoxalin-1-one
MF: $C_9H_5N_3O_2$
FW: 187.2
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 232, 295 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

ODQ is supplied as a crystalline solid. A stock solution may be made by dissolving the ODQ in the solvent of choice, which should be 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 0.5, 15, 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 0.1 mg/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.

Description

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**(2), 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), 1-5 (1996).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/07/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM