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 »

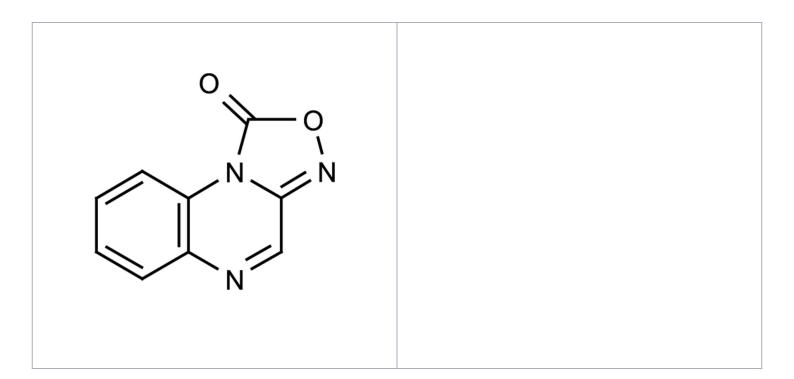
Ordering Information

Order Online »

ALX-270-034-M010	10mg
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

