

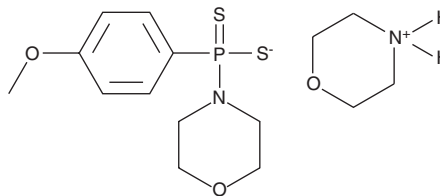
PRODUCT INFORMATION



GY 4137

Item No. 13345

CAS Registry No.: 106740-09-4
Formal Name: (*p*-methoxyphenyl)morpholino-phosphinodithioic acid
MF: C₁₁H₁₆NO₂PS₂ • C₄H₁₀NO
FW: 377.5
Purity: ≥95%
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

GY 4137 is supplied as a crystalline solid. A stock solution may be made by dissolving the GY 4137 in the solvent of choice, which should be purged with an inert gas. GY 4137 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of GY 4137 in these solvents is approximately 15 and 20 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of GY 4137 can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of GY 4137 in PBS (pH 7.2) is approximately 3 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Hydrogen sulfide (H₂S) is a gaseous mediator which, like nitric oxide (NO), has numerous profound actions in mammalian physiology. GY 4137 is a water-soluble, slow-releasing hydrogen sulfide (H₂S) donor.¹ When given intravenously, it demonstrates vasodilator and anti-hypertensive activity in rats, in either the acute (L-NAME-induced) or chronic (spontaneously hypertensive) hypertension models.¹ Intravenous GY 4137 also protects against endotoxic shock in rats, inhibiting tumor necrosis factor- α , interleukin (IL)-1 β , and IL-6 production and reducing NF- κ B activation, iNOS and cyclooxygenase-2 expression, and NO and prostaglandin E₂ generation.²

References

- Li, L., Whiteman, M., Guan, Y.Y., *et al.* Characterization of a novel, water-soluble hydrogen sulfide releasing molecule (GY4137): New insights into the biology of hydrogen sulfide. *Circulation* **117(18)**, 2351-2360 (2008).
- Li, L., Salto-Tellez, M., Tan, C.H., *et al.* GY4137, a novel hydrogen sulfide-releasing molecule, protects against endotoxic shock in the rat. *Free Radic. Biol. Med.* **47(1)**, 103-113 (2009).

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, 03/20/2024

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