PRODUCT INFORMATION



4-hydroxy-3-Methoxyphenylglycol (piperazine salt) Item No. 19714

CAS Registry No.: Formal Name:	67423-45-4 1-(4-hydroxy-3-methoxyphenyl)-		
	1,2-ethanediol compound with piperazine	ОН 	H
Synonyms:	MHPG piperazine, MOPEG piperazine	O OH	• 1/2
MF:	$C_9H_{12}O_4 \bullet_{1/2}C_4H_{10}N_2$		
FW:	227.3	но	N/
Purity:	≥98%		H
UV/Vis.:	λ _{max} : 229, 279 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

4-hydroxy-3-Methoxyphenylglycol (piperazine salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the 4-hydroxy-3-methoxyphenylglycol (piperazine salt) in the solvent of choice, which should be purged with an inert gas. 4-hydroxy-3-Methoxyphenylglycol (piperazine salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 4-hydroxy-3-methoxyphenylglycol (piperazine salt) in these solvents is approximately 10, 15, and 25 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 4-hydroxy-3-methoxyphenylglycol (piperazine salt) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 4-hydroxy-3-methoxyphenylglycol (piperazine salt) in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

4-hydroxy-3-Methoxyphenylglycol (piperazine salt) is a metabolite of norepinephrine derived in part by deamination of normetanephrine or by O-methylation of dihydroxyphenylglycol after cellular uptake of either intermediate.¹ It is released into the blood and cerebrospinal fluid and has been used as a biochemical index of recent sympathetic nervous system activity.²

References

- 1. Kopin, I.J., Axelrod, J., and Gordon, E. The metabolic fate of H3-epinephrine and C14-metanephrine in the rat. J. Biol. Chem. 236, 2109-2136 (1961).
- Goldstein, D. S., Eisenhofer, G., and Kopin, I. J. Sources and significance of plasma levels of catechols and 2. their metabolites in humans. J. Pharm. Exp. Ther. 305(3), 800-811 (2003).

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

SAFFTY 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

uyer 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, 10/20/2022

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