

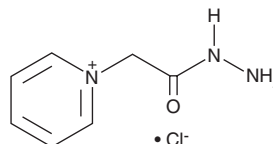
PRODUCT INFORMATION



Girard's Reagent P

Item No. 601541

CAS Registry No.: 1126-58-5
Formal Name: 1-(2-hydrazinyl-2-oxoethyl)-pyridinium, monochloride
Synonyms: Girard P hydrazine, GirP, GP
MF: C₇H₁₀N₃O • Cl
FW: 187.6
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

Girard's reagent P is supplied as a crystalline solid. A stock solution may be made by dissolving the Girard's reagent P in the solvent of choice, which should be purged with an inert gas. Girard's reagent P is soluble in DMSO at a concentration of approximately 1 mg/ml.

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 Girard's reagent P can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of Girard's reagent P in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Girard's reagent P is a cationic hydrazine reagent.¹ It has been used as a derivatization reagent to quantify glycans, oxysterols, and 5-methylcytosine derivatives.²⁻⁴

References

1. Wheeler, O.H. The Girard Reagents. *Chem. Rev.* **62(3)**, 205-221 (1962).
2. Wang, C., Wu, Z., Yuan, J., *et al.* Simplified quantitative glycomics using the stable isotope label Girard's reagent P by electrospray ionization mass spectrometry. *J. Proteome Res.* **13(2)**, 372-384 (2014).
3. Griffiths, W.J., Abdel-Khalik, J., Crick, P.J., *et al.* Sterols and oxysterols in plasma from Smith-Lemli-Opitz syndrome patients. *J. Steroid Biochem. Mol. Biol.* **169**, 77-87 (2017).
4. Tang, Y., Xiong, J., Jiang, H.-P., *et al.* Determination of oxidation products of 5-methylcytosine in plants by chemical derivatization coupled with liquid chromatography/tandem mass spectrometry analysis. *Anal. Chem.* **86(15)**, 7764-7772 (2014).

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, 07/06/2023

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