

PRODUCT INFORMATION

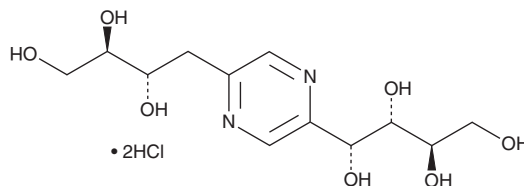


2,5-Deoxyfructosazine (hydrochloride)

Item No. 18662

CAS Registry No.: 17460-13-8
Formal Name: (1R,2S,3R) 1-[5-[(2S,3R)-2,3,4-trihydroxybutyl]-2-pyrazinyl]-1,2,3,4-butanetetrol, dihydrochloride

Synonym: NSC 270912
MF: C₁₂H₂₀N₂O₇ • 2HCl
FW: 377.2
Purity: ≥98%
UV/Vis.: λ_{max}: 274 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

2,5-Deoxyfructosazine (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the 2,5-deoxyfructosazine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. 2,5-Deoxyfructosazine (hydrochloride) is soluble in the organic solvent DMSO. The solubility of 2,5-deoxyfructosazine (hydrochloride) in DMSO is approximately 3 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 2,5-deoxyfructosazine (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 2,5-deoxyfructosazine (hydrochloride) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

2,5-Deoxyfructosazine is a pyrazine derivative that can be found in cured tobacco and is used as a flavoring agent in the food and tobacco industry.^{1,2} Pyrazines, including 2,5-deoxyfructosazine, can be produced either by pyrolytic decomposition of natural compounds or by the reaction of sugars with alcoholic ammonia.² 2,5-Deoxyfructosazine is also formed by the breakdown of D-glucosamine at neutral pH.^{3,4} Like glucosamine, 2,5-deoxyfructosazine has DNA strand breakage activity and strongly inhibits IL-2 production by *Jurkat* cells stimulated with phytohemagglutinin (IC₅₀ = ~1.25 mM).³⁻⁵

References

1. Henry, N., Delépée, R., Seigneuret, J.M., et al. *Talanta* **99**, 816-823 (2012).
2. Rodgman, A. and Perfetti, T.A. *The Chemical Components of Tobacco and Tobacco Smoke*, Second Edition. CRC Press (2013).
3. Zhu, A., Huang, J.-B., Clark, A., et al. *Carbohydr. Res.* **342**(18), 2745-2749 (2007).
4. Shimamura, T., Takamori, A., Ukeda, H., et al. *Biosci. Biotechnol. Biochem.* **67**(2), 295-299 (2003).
5. Sumoto, K., Irie, M., Mibu, N., et al. *Chem. Pharm. Bull. (Tokyo)* **39**(3), 792-794 (1991).

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, 12/15/2022

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