

PRODUCT INFORMATION



D-Allulose-6-phosphate (sodium salt)

Item No. 9004811

Formal Name: D-psicose, 6-(dihydrogen phosphate), monosodium salt

Synonyms: A6P, Allu6P, D-Psicose-6-phosphate

MF: $C_6H_{12}O_9P \cdot Na$

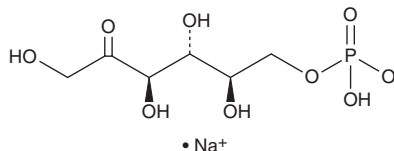
FW: 282.1

Purity: $\geq 90\%$

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

D-Allulose-6-phosphate (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the D-allulose-6-phosphate (sodium salt) in water. D-Allulose-6-phosphate (sodium salt) is slightly soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

D-Allulose-6-phosphate is a monosaccharide. It has been used as a precursor in the enzymatic synthesis of D-allulose, a low-calorie sweetener.¹ D-Allulose-6-phosphate has also been used in the synthesis of type B ribose-5-phosphate isomerase inhibitors.²

References

1. Zhang, Y., Zhou, Z., Luan, H., et al. Advances in the biosynthesis of D-allulose. *World J. Microbiol. Biotechnol.* **40**(12), 375 (2024).
2. Mariano, S., Roos, A.K., Mowbray, S.L., et al. Competitive inhibitors of type B ribose 5-phosphate isomerases: Design, synthesis and kinetic evaluation of new D-allose and D-allulose 6-phosphate derivatives. *Carbohydr. Res.* **344**(7), 869-880 (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, 09/09/2025

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