

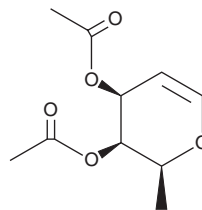
PRODUCT INFORMATION



3,4-Di-O-acetyl-L-fucal

Item No. 30864

CAS Registry No.: 54621-94-2
Formal Name: 2,6-anhydro-1,5-dideoxy-L-arabino-hex-5-enitol, 3,4-diacetate
Synonym: L-Fucal diacetate
MF: C₁₀H₁₄O₅
FW: 214.2
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

3,4-Di-O-acetyl-L-fucal is supplied as a crystalline solid. A stock solution may be made by dissolving the 3,4-di-O-acetyl-L-fucal in the solvent of choice, which should be purged with an inert gas. 3,4-Di-O-acetyl-L-fucal is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 3,4-di-O-acetyl-L-fucal in these solvents is approximately 30 mg/ml.

Description

3,4-Di-O-acetyl-L-fucal is a glycal precursor that has been used in the palladium-catalyzed stereoselective synthesis of glycosides.¹ It has also been used as a precursor in the copper-catalyzed stereoselective synthesis of 2-deoxyglycosides.²

References

1. Pal, K.B., Lee, J., Das, M., *et al.* Palladium(II)-catalyzed stereoselective synthesis of C-glycosides from glycals with diaryliodonium salts. *Org. Biomol. Chem.* **18(12)**, 2242-2251 (2020).
2. Palo-Nieto, C., Sau, A., Jeanneret, R., *et al.* Copper reactivity can be tuned to catalyze the stereoselective synthesis of 2-deoxyglycosides from glycals. *Org. Lett.* **22(5)**, 1991-1996 (2020).

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, 05/13/2020

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