

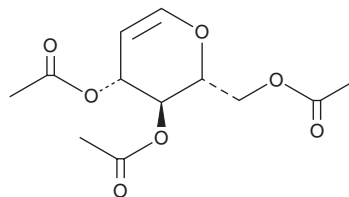
# PRODUCT INFORMATION



## 3,4,6-Tri-O-acetyl-D-glucal

Item No. 39976

**CAS Registry No.:** 2873-29-2  
**Formal Name:** 3,4,6-triacetate 1,5-anhydro-2-deoxy-D-arabino-hex-1-enitol  
**Synonym:** Tri-O-acetyl-D-glucal  
**MF:** C<sub>12</sub>H<sub>16</sub>O<sub>7</sub>  
**FW:** 272.3  
**Purity:** ≥98%  
**Supplied as:** A 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

3,4,6-Tri-O-acetyl-D-glucal is supplied as a solid. A stock solution may be made by dissolving the 3,4,6-tri-O-acetyl-D-glucal in the solvent of choice, which should be purged with an inert gas. 3,4,6-Tri-O-acetyl-D-glucal is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

### Description

3,4,6-Tri-O-acetyl-D-glucal is a derivative of D-(+)-glucose (Item No. 23733) and a precursor in the synthesis of a variety of compounds, including in the synthesis of cyclic ketene acetals and the enantioselective synthesis of spliceostatin A and thailanstatin A derivatives with spliceosome inhibitory activity.<sup>1-3</sup>

### References

1. Hardy, C., Levere, M.E., Kociok-Köhn, G., *et al.* Radical ring opening polymerization of cyclic ketene acetals derived from D-glucal. *ACS Macro Lett.* **12(11)**, 1443-1449 (2023).
2. Ghosh, A.K., Reddy, G.C., Kovala, S., *et al.* Enantioselective synthesis of a cyclopropane derivative of spliceostatin A and evaluation of bioactivity. *Org. Lett.* **20(22)**, 7293-7297 (2018).
3. Ghosh, A.K., Veitschegger, A.M., Nie, S., *et al.* Enantioselective synthesis of thailanstatin A methyl ester and evaluation of *in vitro* splicing inhibition. *J. Org. Chem.* **83(9)**, 5187-5198 (2018).

#### 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, 05/06/2024

#### 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