

# PRODUCT INFORMATION



## Acetylated Carbohydrate Mixture 1

Item No. 29367

**Supplied as:** A solution in chloroform

**Storage:** -20°C

**Stability:** ≥1 year

**Concentration:** 50 mg/ml

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

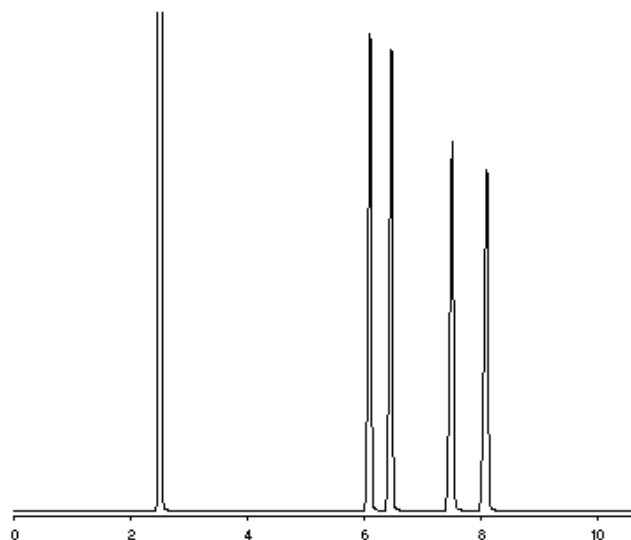
### Description and Data

Acetylated Carbohydrate Mixture 1 contains four pentaacetate alditols. It is intended for use as a reference standard for the detection of plant cell wall polysaccharides, among others, in various sample types by GC-MS. This mixture is supplied in an amber ampule in which the headspace has been purged with nitrogen/argon to prevent lipid oxidation.

Included are:

Elution Order*	Item Name
1	Rhamnitol pentaacetate
2	Fucitol pentaacetate
3	Ribitol pentaacetate
4	Arabinitol pentaacetate

\*GC Column Used: SP-2330, 30 x 0.25 mm x 0.2 μm



#### 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, 10/24/2019

#### 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