

PRODUCT INFORMATION



Plant Sterol Mixture

Item No. 29364

Supplied as: A solution in chloroform

Storage: -20°C

Stability: ≥2 years

Concentration: 25 mg/ml

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

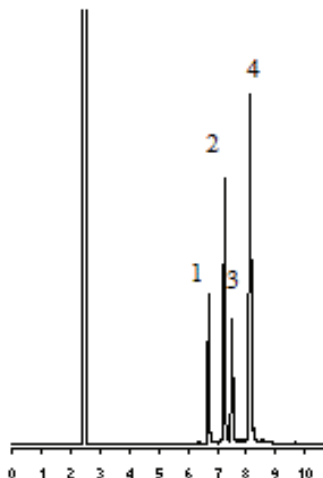
Description and Data

Plant sterols are components of plant cell membranes and occur as free sterols, sterol esters, sterol glycosides, and acylated sterol glycosides.¹ They are precursors of the brassinosteroid family of plant growth regulators and have roles in transmembrane signal transduction and cell proliferation and differentiation. Plant Sterol Mixture contains brassicasterol (14%), campesterol (28%), stigmasterol (12%), and β-sitosterol (43%) and is intended for use as a reference standard for the detection of these plant sterols in various sample types by GC-FID, GC-MS, or LC-MS. This mixture is supplied in an amber ampule in which the headspace has been purged with argon to prevent lipid oxidation.

Included are:

Elution Order*	Item Name
1	Brassicasterol (14%)
2	Campesterol (28%)
3	Stigmasterol (12%)
4	β-Sitosterol (43%)

*GC Column Used: RTX-1, 30 m x 0.25 mm x 0.2 μm



Reference

1. Valitova, J.N., Sulkarnayeva, A.G., and Minibayeva, F.V. Plant sterols: Diversity, biosynthesis, and physiological functions. *Biochemistry (Mosc.)* **81(8)**, 819-834 (2016).

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, 04/28/2026

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