

# PRODUCT INFORMATION



## Plant Sterols Kit

Item No. 28643

**Supplied as:** A solution in chloroform  
**Storage:** -20°C  
**Stability:** ≥2 years

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

### Description

Plant sterols are cell membrane lipids that are found as free sterols, sterol esters, sterol glycosides, and acylsterol glycosides.<sup>1</sup> They are precursors of the brassinosteroid family of plant growth regulators and have roles in transmembrane signal transduction and cell proliferation and differentiation. Plant Sterols Kit contains the following seven individually packaged plant sterols:  $\beta$ -sitosterol, desmosterol, lanosterol, stigmasterol, ergosterol, coprostanol, and cholestanol, as well as three mixtures: steryl glucosides, esterified steryl glucosides, and plant sterol mixture. Steryl glucosides is a mixture of non-esterified  $\beta$ -sitosterol glucoside,  $\beta$ -campesterol glucoside,  $\beta$ -stigmasteryl glucoside, and  $\beta$ - $\Delta$ -5-avenosteryl glucoside. Esterified steryl glucosides contains esterified forms of  $\beta$ -sitosterol glucoside,  $\beta$ -campesterol glucoside, and  $\beta$ -stigmasteryl glucoside. Plant sterol mixture contains brassicasterol (14%), campesterol (28%), stigmasterol (12%), and  $\beta$ -sitosterol (43%). Plant Sterols Kit is intended for use as a reference standard for the detection of these plant sterols in various sample types by TLC, LC-MS, or GC-MS.

### 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, 10/24/2022

#### 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