

CERTIFICATE of ANALYSIS



Plant Sterols Kit

Item No. 28643 • Batch No. 0660803

Purity Specification:

Molecular Formula.:

CAS Number: (unspecified)

Formula Weight: 0

Expiry date: 22SEP2024

Overview

Tests	Results
Egosterol (90%)	25mg
beta-Sitosterol (50%)	100mg
Cholesterol	100mg
Lanosterol (55%)	100mg
Desmosterol	2mg
Coprostanol	5mg
Steryl Glucosides	25mg
Stigmasterol	25mg
Esterified Steryl Glucosides	10mg
*purity 98% unless otherwise stated in brackets	98%
Plant sterol Mixture	25mg/ml in Chloroform

Reviewed and approved by: Justine Albertson

WARNING

THIS PRODUCT IS FOR RESEARCH USE - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE. IT IS THE RESPONSIBILITY OF THE PURCHASER TO DETERMINE SUITABILITY FOR OTHER APPLICATIONS.

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, 10/12/2018

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