

# PRODUCT INFORMATION



## ***Withania somnifera* (Ashwagandha) Aqueous Methanol Extract** Item No. 43251

**Supplied as:** A solution in methanol

**Storage:** -20°C

**Stability:** ≥2 years

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

### **Description**

---

*Withania somnifera* (Ashwagandha) aqueous methanol extract is an extract of the medicinal plant *W. somnifera* containing a crude mixture of bioactive natural products, including withanolide A (Item No. 11797), 12-deoxywithastramonolide (Item No. 11669), and withaferin A (Item No. 11352), among others.

#### **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, 06/04/2025

#### **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](http://WWW.CAYMANCHEM.COM)