

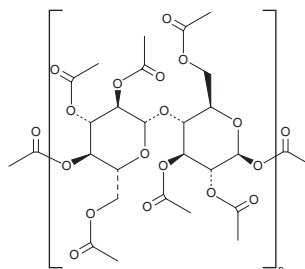
# PRODUCT INFORMATION



## Cellulose triacetate

Item No. 16521

**CAS Registry No.:** 9012-09-3  
**Formal Name:** triacetate cellulose  
**Synonym:** Triacetylcellulose  
**FW:** 678.6  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Description

Cellulose triacetate is a polymer film derived from cellulose treated with an acetylating agent, which renders the cellulose polymers more soluble in organic solvents.<sup>1</sup> Cellulose triacetate can be used to manufacture high-performance membranes characterized by well-defined pores for dialysis and reverse osmosis and is often combined with plasticizer and an ionophore for ion sensing and separation.<sup>1,2</sup> It is also implemented in the production of photographic film and polarized film for liquid crystal displays.<sup>1</sup>

### References

1. Sunohara, T. and Masuda, T. Cellulose triacetate as a high-performance membrane. *Contrib. Nephrol.* **173**, 156-163 (2011).
2. Arous, O., Saad Saoud, F., and Kerdjoudj, H. Cellulose triacetate properties and their effect on the thin films morphology and performance. *IOP Conf. Ser.: Mater. Sci. Eng.* **12**, 012001 (2010).

#### 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, 12/22/2022

#### 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