

PRODUCT INFORMATION



Methylcellulose

Item No. 44711

CAS Registry No.: 9004-67-5
Formal Name: cellulose, methyl ether
Supplied as: A solid
Storage: Room temperature
Stability: ≥4 years

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

Laboratory Procedures

Methylcellulose is supplied as a solid. A stock solution may be made by dissolving the methylcellulose in water. We do not recommend storing the aqueous solution for more than one day.

Description

Methylcellulose is a derivative of cellulose.¹⁻³ Formulations containing methylcellulose have been used as surgical reagents, food additives, adhesives, dispersing agents, emollient laxatives, and as a binder in drug formulations. This product is also available as part of Cayman's Tissue Cryopreservation Reagent Kit (MEDY) (Item No. 502885).

References

1. Nassato, P.L., Pignon, F., Silveira, J.L.M., *et al.* Methylcellulose, a cellulose derivative with original physical properties and extended applications. *Polymers* **7**(5), 777-803 (2015).
2. Searles, G.E. and DesGroseilliers, J.-P. Excipients in topical corticosteroid preparations in Canada. *CMAJ* **141**(5), 396-405 (1989).
3. Lorenzetti, O.J. The influence of epinephrine and pilocarpine upon intraocular pressure following single and combination installation. *Ophthalm. Res.* **2**(6), 328-336 (1971).

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, 11/06/2025

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