

# PRODUCT INFORMATION



## Amylose

Item No. 17664

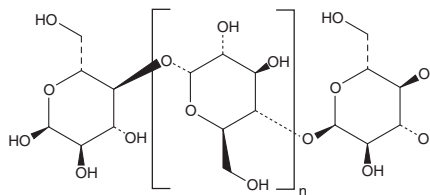
CAS Registry No.: 9005-82-7

MF:  $(C_6H_{10}O_5)_n$

Supplied as: A solid

Storage:  $-20^{\circ}C$

Stability:  $\geq 4$  years



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

### Laboratory Procedures

Amylose is supplied as a solid. A stock solution may be made by dissolving the amylose in the solvent of choice. Amylose is soluble in 0.5 M sodium hydroxide at a concentration of approximately 1 mg/ml.

Amylose is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

### Description

Amylose is a polysaccharide made of  $\alpha$ -D-glucose units, bound to each other through  $\alpha(1\rightarrow4)$  glycosidic bonds. It constitutes about 20% of starch, the major storage carbohydrate in plants. It can act as an indicator, turning blue in the presence of iodine, and has also been coupled to magnetic beads for the isolation and purification of maltose-binding protein fusion proteins.<sup>1,2</sup>

### References

1. Ovecka, M., Bahaji, A., Muñoz, F.J., *et al.* A sensitive method for confocal fluorescence microscopic visualization of starch granules in iodine stained samples. *Plant Signal. Behav.* **7(9)**, 1146-1150 (2012).
2. Affinity chromatography: Methods and protocols, *in* Methods in molecular biology. Zachariou, M., editor, Humana Press, Vol. 421, Second Ed. (2008).

#### 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, 10/19/2022

#### 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