

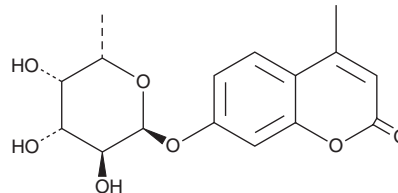
# PRODUCT INFORMATION



## 4-Methylumbelliferyl- $\alpha$ -L-Fucopyranoside

Item No. 39552

**CAS Registry No.:** 54322-38-2  
**Formal Name:** 7-[[6-deoxy- $\alpha$ -L-galactopyranosyl]oxy]-4-methyl-2H-1-benzopyran-2-one  
**Synonyms:** 4-Methylumbelliferyl- $\alpha$ -L-Fucoside, 4-MU- $\alpha$ -L-Fucopyranoside  
**MF:** C<sub>16</sub>H<sub>18</sub>O<sub>7</sub>  
**FW:** 322.3  
**Purity:**  $\geq$ 98%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years



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

### Laboratory Procedures

4-Methylumbelliferyl- $\alpha$ -L-fucopyranoside is supplied as a solid. A stock solution may be made by dissolving the 4-methylumbelliferyl- $\alpha$ -L-fucopyranoside in the solvent of choice, which should be purged with an inert gas. 4-Methylumbelliferyl- $\alpha$ -L-fucopyranoside is soluble in the organic solvent dimethyl formamide.

### Description

4-Methylumbelliferyl- $\alpha$ -L-fucopyranoside is a fluorogenic substrate for  $\alpha$ -L-fucosidase.<sup>1</sup> Upon enzymatic cleavage by  $\alpha$ -L-fucosidase, 4-methylumbelliferone (4-MU) is released and its fluorescence can be used to quantify  $\alpha$ -L-fucosidase activity. 4-MU displays excitation/emission maxima of 350/440 nm, respectively.

### Reference

1. Robinson, D. and Thorpe, R. Fluorescent assay of  $\alpha$ -L-fucosidase. *Clin. Chim. Acta* **55(1)**, 65-69 (1974).

#### 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, 10/16/2023

#### 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