

## 4-Methylumbelliferyl alpha-D-galactopyranoside

Catalog Number: 14017

Unit Size: 10 mg

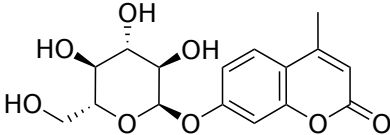
### Product Details

---

Storage Conditions	N/A
Expiration Date	12 months upon receiving

### Chemical Properties

---

Appearance	Solid white
Molecular Weight	338.31
Soluble In	DMSO
Chemical Structure	

### Spectral Properties

---

Excitation Wavelength	N/A
Emission Wavelength	N/A

### Applications

---

4-Methylumbelliferyl  $\alpha$ -D-glucopyranoside is used as a fluorogenic substrate for the identification, characterization and kinetic analysis of  $\alpha$ -D-glucosidases. 4-Methylumbelliferyl  $\alpha$ -D-glucopyranoside is used for the high throughput screening of  $\alpha$ -D-glucosidase inhibitors. It is also suitable for the diagnosis of Pompe's disease (also known as acid-maltase disease and glycogen storage disease II) in kidney and leucocytes. Pompe disease is a rare genetic disorder that causes progressive weakness to the heart and skeletal muscles.