

4-Fluororesorcinol *CAS#: 103068-41-3*

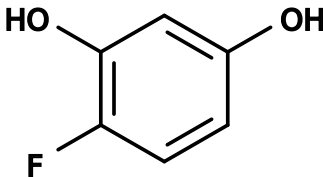
Catalog Number: 4022

Unit Size: 1 g

Product Details

Storage Conditions	Freeze (< -15 °C)
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Solid
Molecular Weight	128.10
Soluble In	DMF
Chemical Structure	

Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

4-Fluororesorcinol is a precursor that can be used for preparing fluorescent dyes such as fluorescein, coumarin and resorufin analogs. It is the key precursor for preparing Oregon Green dyes.