

**2,4-Difluororesorcinol \*CAS 195136-71-1\***

Catalog Number: 4023

Unit Size: 1 g

**Product Details**

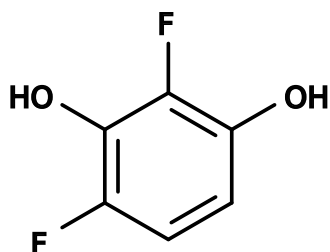
---

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

**Chemical Properties**

---

Appearance	Solid off-white
Molecular Weight	146.09
Soluble In	DMF
Chemical Structure	

**Spectral Properties**

---

Excitation Wavelength	N/A
Emission Wavelength	N/A

**Applications**

---

2,4-Difluororesorcinol 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 Pacific Blue dyes.