

## 3,6-Dichlorobenzene-1,2,4-tricarboxylic anhydride [3,6-Dichlorotrimellitic anhydride]

Catalog Number: 4021

Unit Size: 1 g

### Product Details

---

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

### Chemical Properties

---

Appearance	Solid light gray
Molecular Weight	261.01
Soluble In	DMF
Chemical Structure	

### Spectral Properties

---

Excitation Wavelength	N/A
Emission Wavelength	N/A

### Applications

---

3,6-Dichlorotrimellitic anhydride is the key precursor that is used for preparing a variety of dichlorinated fluoresceins and rhodamines such as TET and HEX. These chlorinated fluoresceins and rhodamines are widely used for labeling oligos and in DNA sequencing.