

Product Information Sheet

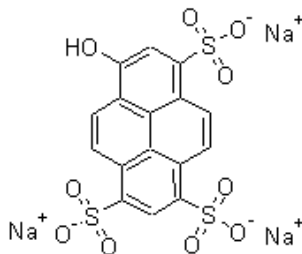
Ordering Information

Product Number: 85
Product Name: 8-Hydroxypyrene-1,3,6-trisulfonic acid, trisodium salt [HPTS]*
Unit Size: 1 g
Storage Conditions: <-15 °C and kept from light and moisture
Expiration Date: 12 months upon receiving

Chemical and Spectral Properties

Molecular Weight: 524.38

Chemical Structure:



Soluble in: Water
Excitation Wavelength: 454 nm
Emission Wavelength: 511 nm

Application Notes

8-Hydroxypyrene-1,3,6-trisulfonic acid, trisodium salt [HPTS] is the cell-impermeable fluorescent pH indicator used for slightly acidic pH range (pKa = 6.4). It is highly water-soluble.