

Product Information Sheet

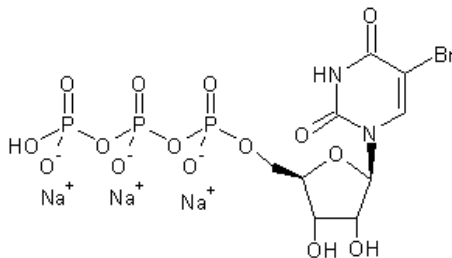
Ordering Information

Product Number: 17032
Product Name: BrUTP [5-Bromouridine 5'-triphosphate] *CAS# 161848-60-8*
Unit Size: 100 µL
Storage Conditions: <-15 °C and kept from light and moisture
Expiration Date: 12 months upon receiving

Chemical and Spectral Properties

Molecular Weight: 628.98

Chemical Structure:



Soluble in: Water

Excitation Wavelength: N/A

Emission Wavelength: N/A

Application Notes

BrUTP can be used for incorporating into RNA for the subsequent detection with anti-BrdU antibodies. BrUTP incorporation into RNA is also a tool for random-mutation introduction.