

Product Information Sheet

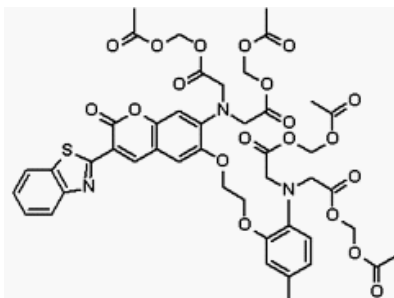
Ordering Information

Product Number: 21054
Product Name: BTC, AM
Unit Size: 1 mg
Storage Conditions: <-15 °C and kept from light and moisture
Expiration Date: 12 months upon receiving

Chemical and Spectral Properties

Molecular Weight: 979.92

Chemical Structure:



Soluble in: DMSO
Excitation Wavelength: 401 nm
Emission Wavelength: 529 nm

Application Notes

This cell-permeant coumarin Ca^{2+} indicator BTC, AM exhibits a shift in excitation maximum from about 480 nm to 400 nm upon binding Ca^{2+} , enabling ratiometric calcium measurements. Due to its high selectivity and low affinity for Ca^{2+} ($K_d \sim 7 \mu\text{M}$) BTC is often used for the quantitation of high intracellular Ca^{2+} levels. In addition, BTC, AM has also been used for monitoring potassium channel since thallium ion enhances the fluorescence of BTC.