

Product Information Sheet

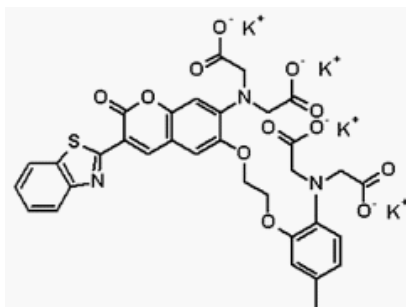
Ordering Information

Product Number: 21053
Product Name: BTC, tetrapotassium salt
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: 844.02

Chemical Structure:



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

Application Notes

This cell-impermeant coumarin Ca^{2+} indicator 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+} ($K_d \sim 7 \text{ uM}$) BTC is often used for the quantitation of high Ca^{2+} levels.