
Product Information Sheet

Ordering Information

Product Number: 20648

Product Name: Calbryte™-520XL-Dextran

Unit Size: 1 mg

Storage Conditions: Freeze (<-15 °C), Desiccated, Avoid Light

Expiration Date: 12 months upon receiving

Chemical and Spectral Properties

Appearance:

Molecular Weight: ~11000

Soluble In: Water

Excitation Wavelength: 492

Emission Wavelength: 514

Application Notes

Calbryte 520XL dextran is a new fluorescent and cell-impermeable calcium indicator with extremely low affinity. It has $K_d \sim 500 \mu\text{M}$ which is similar to the well-known Rhod 5N, but much more stable than Rhod-5N. Calbryte 520XL produces bright fluorescence signal in the presence of calcium ion in high concentration. It has the identical excitation and emission wavelength as Fluo-4, thus the same Fluo-4 assay settings can be readily applied to Calbryte 520XL-based calcium assays. Calbryte 520XL is an excellent alternative to Rhod-5N. We also offer Calbryte 520XL, AM (#20646), potassium salt (#20645), and Calbryte 520XL azide (#20643) that can be readily conjugated to a carrier through the well-know click chemistry.