
Product Information Sheet

Ordering Information

Product Number:	20706
Product Name:	Calbryte™-590, potassium salt
Unit Size:	5x50 ug
Storage Conditions:	Freeze (<-15 °C), Desiccated, Avoid Light
Expiration Date:	12 months upon receiving

Chemical and Spectral Properties

Appearance:	Solid
Molecular Weight:	1074.98
Soluble In:	Water
Excitation Wavelength:	573
Emission Wavelength:	588

Application Notes

The intracellular calcium flux assay is a widely used method in monitoring signal transduction pathways and high throughput screening of G protein-coupled receptors (GPCRs) and calcium channel targets. Followed by Rhod-2 being introduced in 1989, Rhod-4 and Cal-590 were later developed with improved signal/background ratio, and they became the widely used red fluorescent Ca²⁺ indicators for confocal microscopy, flow cytometry and high throughput screening applications. Calbryte 590 is a new generation of red fluorescent indicators for the measurement of intracellular calcium. Its greatly improved signal/background ratio and intracellular retention properties make CalBryte 590 the most robust red fluorescent indicator for evaluating GPCR and calcium channel targets as well as for screening their agonists and antagonists in live cells.