

## Calcein Deep Red™

Catalog Number: 21902

Unit Size: 1 mg

### Product Details

---

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

### Chemical Properties

---

Appearance	Solid black
Molecular Weight	373.66
Soluble In	DMSO

### Spectral Properties

---

Excitation Wavelength	643 nm
Emission Wavelength	663 nm

### Applications

---

Calcein AM is one the most popular fluorescent probes used for labeling and monitoring cellular functions of live cells. However, the single color of Calcein AM makes it impossible to use this valuable reagent in the multicolor applications. For example, it is impossible to use Calcein AM in GFP-transfected cells due to the same color to GFP. To address this color limitation of Calcein AM, we have developed Calcein Orange™, Calcein Red™ and Calcein Deep Red™. These new Calcein AM analogs enable the multicolor labeling and functional analysis of live cells in combination with Calcein AM. Calcein Deep Red™ is the esterase-hydrolyzed product of Calcein Deep Red™ acetate in cells. Calcein Deep Red™ dye can be monitored with the common Cy5 filter set. It is used as a fluorescence reference standard for Calcein Deep Red™ acetate.