

**CytoCalcein™ Violet 660, sodium salt \*Excited at 405 nm\***

Catalog Number: 21903

Unit Size: 1 mg

**Product Details**

---

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

**Chemical Properties**

---

Appearance	Solid
Molecular Weight	662.63
Soluble In	DMSO

**Spectral Properties**

---

Excitation Wavelength	402 nm
Emission Wavelength	660 nm

**Applications**

---

CytoCalcein™ Violet 660 is the esterase-cleaved product of CytoCalcein™ Violet 660 AM in live cells. CytoCalcein™ Violet 660 is not cell-permeable. It is the reference product for the CytoCalcein™ Violet 660 AM that is designed for labeling live cells in the same way as Calcein, AM does. CytoCalcein™ Violet 660 is well excited with the Violet laser at 405 nm, which is now equipped in most flow cytometers. It is also well-excited by the excitation sources of fluorescence microscopes. This exceptional spectral separation from the typical existing FACS fluorophores provides additional options for multiplexing experiments with a flow cytometer or a fluorescence microscope. CytoCalcein™ Violet 660 has been developed for multiplexing flow cytometric and fluorescence imaging applications. It provides a new unique color for flow cytometric analysis of live cells.