

**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.