

## mFluor™ Blue 660 tyramide

Catalog Number: 11051

Unit Size: 200 Slides

### Product Details

---

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

### Chemical Properties

---

|                  |        |
|------------------|--------|
| Appearance       | Solid  |
| Molecular Weight | 770.00 |
| Soluble In       | DMSO   |

### Spectral Properties

---

|                       |        |
|-----------------------|--------|
| Excitation Wavelength | 481 nm |
| Emission Wavelength   | 663 nm |

### Applications

---

mFluor™ Blue 660 tyramide is a fluorescent labeling reagent used in immunofluorescence staining and in situ hybridization. Tyramide is a small molecule that can diffuse through tissue sections or cell membranes and subsequently be enzymatically amplified to produce a localized, highly fluorescent signal. The mFluor™ Blue 660 dye is conjugated to tyramide to create a fluorescently labeled tyramide that can be used for visualizing specific target molecules or structures within biological samples. mFluor™ Blue 660 tyramide has the largest Stokes Shift among all the commercial tyramide reagents. The high sensitivity and specificity of mFluor™ Blue 660 tyramide makes it an excellent choice for fluorescence imaging and detection in various histochemical fluorescence imaging applications.