

**mFluor™ Blue 585 acid**

Catalog Number: 72101

Unit Size: 1 mg

**Product Details**

---

Storage Conditions Freeze (&lt; -15 °C), Minimize light exposure

Expiration Date 24 months upon receiving

**Chemical Properties**

---

Appearance Solid

Molecular Weight 580.64

Soluble In DMSO

**Spectral Properties**

---

Excitation Wavelength 491 nm

Emission Wavelength 578 nm

**Applications**

---

mFluor™ Blue 585 acid is a small molecule fluorescent dye with spectral properties comparable to R-phycoerythrin (RPE)-type conjugates when used with 488 nm excitation and red fluorescence detection. The dye is water-soluble and, when converted into appropriate reactive forms for conjugation, can be used to generate protein conjugates that are efficiently excited by 488 nm laser sources and detected using TRITC-compatible settings. These properties make mFluor™ Blue 585 derivatives suitable for applications such as flow cytometry and fluorescence-based assays.

Compared to protein-based fluorophores such as RPE, mFluor™ Blue 585 dyes are small organic molecules, which can facilitate more efficient conjugation to proteins. As part of AAT Bioquest's mFluor™ dye family, they are designed with large Stokes shifts and are optimized for protein labeling, particularly antibodies. These dyes are characterized by bright fluorescence, good photostability, and minimal quenching on proteins, and have been applied in multicolor workflows, including spectral flow cytometry using common excitation sources (e.g., 350, 405, 488, 555, 633, and 647 nm).