

**iFluor™ 690 succinimidyl ester**

Catalog number: 1042

Unit size: 1 mg

**Product Details**

---

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

**Chemical Properties**

---

Appearance	Solid
Soluble In	DMSO

**Applications**

---

AAT Bioquest's iFluor™ dyes are optimized for labeling proteins, in particular, antibodies. These dyes are bright, photostable and have minimal quenching on proteins. They can be well excited by the major laser lines of fluorescence instruments (e.g., 350, 405, 488, 555, 633 and 647 nm) or other excitation sources. iFluor™ 690 dyes have fluorescence excitation and emission maxima of ~690 nm and ~705 nm respectively. These spectral characteristics make them an excellent alternative to Alexa Fluor® 700 labeling dye (Alexa Fluor® is the trademark of Invitrogen). iFluor™ 690 SE is reasonably stable and shows good reactivity and selectivity with protein amino groups. Our preliminary tests indicated that it has superior FRET when conjugated to PE, making it an excellent acceptor dyes for preparing tandems with APC, PE and PerCP. iFluor™ 690-based tandems provide unique colors that are extremely useful for spectral flow cytometry applications.