

## ReadiCleave™ iFluor® 594 AML-NHS ester

Catalog Number: 7007G2

Unit Size: 1 mg

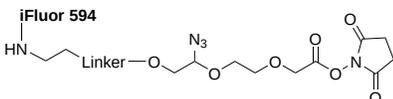
### Product Details

---

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

### Chemical Properties

---

Appearance	Solid purple
Molecular Weight	1622.93
Soluble In	DMSO
Chemical Structure	

### Spectral Properties

---

Excitation Wavelength	587 nm
Emission Wavelength	603 nm

### Applications

---

Many biological molecules can be readily labeled with a fluorescent tag for fluorescence imaging and flow cytometry analysis. However, almost all the existing fluorescent tags are used to permanently labeling biological targets from which the added fluorescent tags cannot be cleaved for further downstream analysis, such as mass spectral analysis or another detection mode. AAT Bioquest's ReadiCleave™ linkers enable fluorescent tags conjugated to a biological target from which the added fluorescent tag can be removed when needed. ReadiCleave™ iFluor® 594 AML uses an azidomethyl linker that can be cleaved with TCEP to remove the iFluor® 594 fluorophore from the target molecule. The cleavage can be carried out by adding 10-100 mM TCEP solution (pH 7.5) and incubating at 65 °C for 1-5 min. iFluor® 594 is a superior replacement to Alexa Fluor® 594. iFluor® 594 and Alexa Fluor® 594 have very similar spectral properties.