

## ReadiCleave™ XFD532 AML-NHS ester

Catalog Number: 7001G2

Unit Size: 1 mg

### Product Details

---

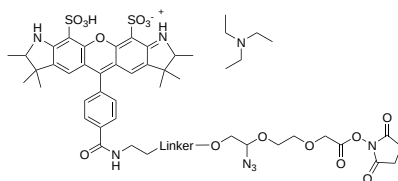
Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

### Chemical Properties

---

Appearance	Solid orange-red
Molecular Weight	1287.47
Soluble In	DMSO

Chemical Structure



### Spectral Properties

---

Excitation Wavelength	534 nm
Emission Wavelength	553 nm

### Applications

---

Fluorescence-based methods have many advantages for biological detections in terms of sensitivity and convenience. Many biological molecules can be readily labeled with a fluorescent tag for fluorescence imaging and flow cytometry analysis. However, most of 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. AAT Bioquest's ReadiCleave™ linkers enable fluorescent tags conjugated to a biological target from which the added fluorescent tag can be removed when needed. This ReadiCleave™ AML XFD532 (mw = 1287.47) contains an azidomethyl linker that can be cleaved with TCEP to remove the XFD532 fluorophore from the target molecule. The cleavage can be carried out by adding 10 mM TCEP solution (pH 7.5), and incubating at 65 °C for 1-5 min. XFD532 is the same fluorophore to Alexa Fluor® 532 (Alexa Fluor® is the trademark of ThermoFisher).