

ReadiCleave™ FITC AML-NHS ester

Catalog Number: 7004G2

Unit Size: 1 mg

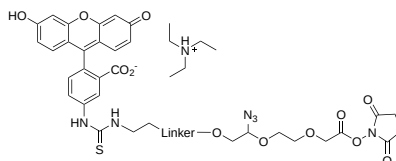
Product Details

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

Chemical Properties

Appearance	Solid yellow
Molecular Weight	1080.14
Soluble In	DMSO

Chemical Structure



Spectral Properties

Excitation Wavelength	498 nm
Emission Wavelength	517 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 FITC (mw = 1080.14) contains an azidomethyl linker that can be cleaved with TCEP to remove the fluorescein 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.