

ReadiCleave[™] iFluor 488 AML-NHS ester

Catalog number: 7005 Unit size: 1 mg

Product Details	
Storage Conditions	Freeze (<-15 °C), Minimize light exposure
Expiration Date	12 months upon receiving
Chemical Properties	
Appearance	Solid
Molecular Weight	1296.43
Soluble In	DMSO
Spectral Properties	
Excitation Wavelength	491 nm
Emission Wavelength	516 nm
A H H	

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 tags conjugated to a biological target from which the added fluorescent tags conjugated to a biological target from which the added fluorescent tag can be removed when needed. This ReadiCleave™ AML iFluor™ 488 contains an azidomethyl linker that can be cleaved with TCEP to remove the iFluor™ 488 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. iFluor™ 488 is a superior replacement to Alexa Fluor® 488. iFluor™ 488 and Alexa Fluor® 488 have very similar spectral properties.