

ReadiCleave™ AML Cy5 NHS ester

Catalog Number: 7000G2

Unit Size: 1 mg

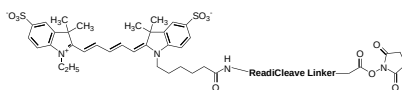
Product Details

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

Chemical Properties

Appearance	Solid blue
Molecular Weight	1317.58
Soluble In	DMSO

Chemical Structure



Spectral Properties

Excitation Wavelength	651 nm
Emission Wavelength	670 nm

Applications

Fluorescence-based methods have many advantages for biological detections in terms of sensitivity. Many biological molecules are labeled with fluorescent tags 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. 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 Cy5 (mw = 1317.58) contains an azidomethyl linker that can be cleaved with TCEP to remove the Cy5 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.