

ReadiCleave™ iFluor 700 SSL-NHS ester

Catalog number: 7053 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 1140.41

Soluble In DMSO

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 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™ 700 succinimidyl ester contains a S-S linker that can be cleaved with TCEP or DTT to remove the iFluor™ 700 fluorophore from the target molecule. iFluor™ 700 has almost identical spectra of Alexa Fluor 700 with higher fluorescence quantum yield and improved stability. The cleavage can be carried out by adding DTT or TCEP solution and incubating from room temperature to 65 °C for 1-10 min.