

Annexin V-iFluor® 568 conjugate

Catalog Number: 20078

Unit Size: 100 tests

Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	6 months upon receiving

Chemical Properties

Appearance	Liquid
Molecular Weight	N/A
Soluble In	Water

Spectral Properties

Excitation Wavelength	568 nm
Emission Wavelength	587 nm

Applications

Annexin V-iFluor® 568 Conjugate is a fluorescent apoptosis detection reagent that combines Annexin V with iFluor® 568, a bright orange-red fluorophore with excitation/emission maxima of ~566/586 nm. This conjugate enables sensitive and direct detection of phosphatidylserine (PS) exposure on the outer plasma membrane—a hallmark of early apoptosis.

By binding PS in a calcium-dependent manner, Annexin V-iFluor® 568 serves as a robust tool for identifying apoptotic cells before morphological changes occur. It is suitable for fluorescence microscopy, flow cytometry, and high-content screening of cell death pathways across a range of research applications.