

ReadiLink™ Rapid XFD750 Antibody Labeling Kit *XFD750 Same Structure to Alexa Fluor™ 750*Catalog number: 1279
Unit size: 2x50 ug Labelings**Product Details**

Expiration Date 6 months upon receiving

Spectral Properties

Excitation Wavelength 752 nm

Emission Wavelength 776 nm

Applications

XFD750 is manufactured by AAT Bioquest, and it has the same chemical structure of Alexa Fluor® 750 (Alexa Fluor® is the trademark of ThermoFisher). Readilink™ protein labeling technology is the most robust and convenient tool for preparing fluorescent antibody conjugates used for fluorescence imaging and flow cytometry applications. ReadiLink™ Rapid XFD750 Antibody Labeling Kit provides the most convenient tool for making XFD750-labeled antibody conjugates. XFD750 dye used in this ReadiLink™ kit is reasonably stable and shows good reactivity and selectivity with protein amino groups. The kit has all the essential components for labeling ~2x50 ug antibody. Each of the two vials of XFD750 dye provided in the kit is optimized for labeling ~50 µg antibody. ReadiLink™ Rapid XFD750 Antibody Labeling Kit requires minimal hands-on time. The prepared XFD750 conjugates (with the kit) are ready to use for fluorescence imaging and flow cytometry applications without further purifications needed. It provides the most convenient method to prepare XFD750-labeled antibodies.