

**ReadiLink™ xtra Rapid XFD555 Antibody
Labeling Kit *BSA-Compatible, XFD555 Same
Structure to Alexa Fluor™ 555***Catalog number: 1980
Unit size: 2 Labelings**Product Details**

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

Spectral Properties

Excitation Wavelength	553 nm
Emission Wavelength	568 nm

Applications

XFD555 is manufactured by AAT Bioquest, and it has the same chemical structure of Alexa Fluor® 555 (Alexa Fluor® is the trademark of ThermoFisher). ReadiLink™ xtra rapid antibody labeling kits require essentially only 2 simple mixing steps without a column purification needed. Specially formulated and preactivated XFD555 (chemically equivalent to Alexa Fluor® 555) used in this ReadiLink™ kit is quite stable and shows good reactivity and selectivity with antibodies. The kit has all the essential components for labeling ~2x50 ug antibody. Each of the two vials of preactivated XFD555 dye provided in the kit is optimized for labeling ~50 µg antibody. ReadiLink™ xtra XFD555 rapid antibody labeling kit provides a convenient and robust method to label monoclonal and polyclonal antibodies with red fluorescent XFD555 fluorophore. XFD555 is one of the most used fluorophores for labeling antibodies.