

## ReadiLink™ Rapid XFD647 Antibody Labeling Kit

Catalog Number: 1278

Unit Size: 2x50 ug Labelings

### Product Details

---

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
--------------------	--

Expiration Date	6 months upon receiving
-----------------	-------------------------

### Chemical Properties

---

Appearance	Solid
------------	-------

Molecular Weight	N/A
------------------	-----

Soluble In	DMSO
------------	------

### Spectral Properties

---

Excitation Wavelength	650 nm
-----------------------	--------

Emission Wavelength	671 nm
---------------------	--------

### Applications

---

XFD647 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 647 (Alexa Fluor® is the trademark of Thermo Fisher). 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 XFD647 Antibody Labeling Kit provides the most convenient tool for making XFD647-labeled antibody conjugates. XFD647 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 XFD647 dye provided in the kit is optimized for labeling ~50 µg antibody. ReadiLink™ Rapid XFD647 Antibody Labeling Kit requires minimal hands-on time. The prepared XFD647 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 XFD647-labeled antibodies.