

Wheat Germ Agglutinin, XFD594 Labeled *XFD594 Same Structure to Alexa Fluor™ 594*

Catalog number: 25509 Unit size: 1 mg

Product Details	
Storage Conditions	Freeze (<-15 °C), Minimize light exposure
Expiration Date	12 months upon receiving
Chemical Properties	
Appearance	Solid
Soluble In	Water
Spectral Properties	
Excitation Wavelength	590 nm
Emission Wavelength	618 nm
Applications	

XFD594 is manufactured by AAT Bioquest, and it has the same chemical structure of Alexa Fluor® 594 (Alexa Fluor® is the trademark of ThermoFisher). Wheat germ agglutinin (WGA) is a lectin that binds to N-acetyl-D-glucosamine and sialic acid. Since WGA binds to glycoconjugates its derivatives and conjugates are widely used to label cell membranes and fibrotic scar tissue for fluorescence imaging and analysis. XFD594 conjugate of WGA is equivalent to the Alexa Fluor® 594 conjugate of WGA. It exhibits the bright, red fluorescence. XFD594 WGA conjugate binds to sialic acid and N-acetylglucosaminyl residues.