

Wheat Germ Agglutinin, XFD594 Labeled

Catalog Number: 25509

Unit Size: 1 mg

Product Details

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

Expiration Date	12 months upon receiving
-----------------	--------------------------

Unit Details

Units	25509 (1 mg)
-------	--------------

Reconstitution Volume	0.5 mL ddH ₂ O
-----------------------	---------------------------

Chemical Properties

Appearance	Solid purple
------------	--------------

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

Soluble In	Water
------------	-------

Spectral Properties

Excitation Wavelength	590 nm
-----------------------	--------

Emission Wavelength	618 nm
---------------------	--------

Applications

XFD594 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 594 (Alexa Fluor® is the trademark of Thermo Fisher). 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 bright, red fluorescence. XFD594 WGA conjugate binds to sialic acid and N-acetylglucosaminyl residues.