

Wheat Germ Agglutinin, XFD532 Labeled

Catalog Number: 25502

Unit Size: 1 mg

Product Details

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

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

Unit Details

Units	25502 (1 mg)
-------	--------------

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

Chemical Properties

Appearance	Solid pink
------------	------------

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

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

Spectral Properties

Excitation Wavelength	534 nm
-----------------------	--------

Emission Wavelength	553 nm
---------------------	--------

Applications

XFD532, manufactured by AAT Bioquest, has a similar chemical structure to Alexa Fluor® 532, a trademark of Thermo Fisher. Wheat germ agglutinin (WGA) is a well-researched lectin known for its valuable biological applications. Due to its ability to bind to glycoconjugates, WGA derivatives, and conjugates are widely used to label cell membranes and fibrotic scar tissue for fluorescence imaging and analysis. WGA specifically targets sequences of β -1,4-GlcNAc-linked residues known as chitodextrins. Each monomer possesses two identical, non-interacting binding sites that complement 3 or 4 β -1,4-GlcNAc units. Among the tested monosaccharides, only GlcNAc exhibits binding to WGA, while ManNAc does not bind, and GalNAc demonstrates weak binding. XFD532 WGA conjugate, like its counterpart Alexa Fluor® 532 WGA conjugate, emits a bright yellow fluorescence and is useful in a variety of applications, including immunofluorescence (IF), immunohistochemistry (IHC), and flow cytometry (FC).