

Wheat Germ Agglutinin, XFD555 Labeled

Catalog Number: 25503

Unit Size: 1 mg

Product Details

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

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

Unit Details

Units	25503 (1 mg)
-------	--------------

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

Chemical Properties

Appearance	Solid red
------------	-----------

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

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

Spectral Properties

Excitation Wavelength	553 nm
-----------------------	--------

Emission Wavelength	568 nm
---------------------	--------

Applications

XFD555, manufactured by AAT Bioquest, has a similar chemical structure to Alexa Fluor® 555, 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 yeast bud scars, fibrotic scar tissue, and the cell membrane of gram bacteria and mammalian cells. 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. XFD555 WGA conjugate, like its counterpart Alexa Fluor® 555 WGA conjugate, emits a bright orange fluorescence and is useful in a variety of applications, including immunofluorescence (IF), immunohistochemistry (IHC), and flow cytometry (FC).