

Product Information

Andy Fluor™ 488 WGA

Catalog Number	Unit Size
C015	5 mg

Storage upon receipt:

- -20°C
- Protect from light

Ex/Em: 495/520 nm

Product Description

Wheat germ agglutinin (WGA) is one of the most widely used lectins in cell biology. Our Andy Fluor™ 488 WGA conjugate is useful probe for detecting glycoconjugates, i.e., N-acetylglucosamine and N-acetylneuraminic acid (sialic acid) residues, on cell membranes. Andy Fluor™ 488 WGA can be used in fixed and live cells that have not been permeabilized, to visualize cell membranes. Andy Fluor™ 488 WGA is also used as a retrograde tracer for neuronal mapping where it will also cross synapses.

Application

Cell membrane stain.