

Wheat Germ Agglutinin (WGA) Conjugates

A wheat germ agglutinin (WGA) lectin labeled with Biotium's CF® Dyes, biotin, or HRP. It is a carbohydrate-binding lectin that has high affinity for sialic acid and N-acetylglucosamine and stains yeast bud scars as well as the cell membrane of gram+ bacteria and mammalian cells.



Product Description

Wheat germ agglutinin (WGA) is a carbohydrate-binding lectin that has high affinity for sialic acid and N-acetylglucosamine moieties of glycoproteins. As a result, WGA conjugates will label yeast bud scars and the cell membrane of gram bacteria and mammalian cells.

- Stain yeast bud scars as well as the cell membrane of gram bacteria and mammalian cells or tissues
- Suitable for Western Blot, ELISA, immunohistochemistry and other standard immunoassay applications
- A retrograde or anterograde neuronal tracer
- Withstands fixation and permeabilization
- Choice of 13 CF® Dyes from UV to near-infrared, biotin, or HRP
- Superior CF® Dyes are bright, photostable, and water-soluble

WGA is commonly used to label glycoproteins for imaging of the plasma membrane in live or fixed cells, for staining of tissue sections, or for western blotting. WGA can be used as a gram stain to fluorescently label gram bacteria but not gram- bacteria. WGA also binds to the bud scars on budding yeast such as *Saccharomyces cerevisiae*.

Find the Right Stain for Your Application

WGA and other lectins are carbohydrate binding proteins that recognize specific sugar moieties on glycoproteins. The presence and distribution of these targets vary between cell types and tissues. As a result, other [cell surface stains](#) or other lectin conjugates, [Concanavalin A \(Con A\) Conjugates](#) and [PNA Lectin Conjugates](#), may produce better surface staining and may be more appropriate for your cell type. Lectin conjugates can be used to selectively stain the cell surface of live cells, and withstand fixation and permeabilization. When cells are fixed and permeabilized before staining, fluorescent lectins stain both cell surface and organelles in the secretory pathway. Lectins may be toxic or stimulatory to live cells depending on cell type. To find the right stain for your application, see our [Membrane & Cell Surface Stains Comparison](#). See our [Cellular Stains Table](#) for more information on how our dyes stain various organisms.

Superior CF® Dyes

Biotium's next-generation CF® Dyes were designed to be highly water-soluble with advantages in brightness and photostability compared to Alexa Fluor®, DyLight®, and other fluorescent dyes. Learn more about [CF® Dyes](#).

Note: Conjugates of blue-fluorescent dyes like CF®350, CF®405S and CF®405M are not recommended for detecting low abundance targets and may be challenging to use in tissue specimens. Blue dyes have lower fluorescence and photostability, and cells and tissue have high autofluorescence in blue wavelengths, resulting in lower signal to noise compared to other colors.

Wheat Germ Agglutinin Conjugates

Full List of Lectin Conjugates

CF is a registered trademark of Biotium, Inc. Alexa Fluor, Texas Red, and DyLight are registered trademarks of Thermo Fisher Scientific.

References

Download a list of curated [CF® Dye references](#).

Product attributes

Probe cellular localization	Membrane/cell surface
For live or fixed cells	For fixed cells, For live/intact cells
Cell permeability	Membrane impermeant
Fixation options	Fix before staining (formaldehyde), Fix after staining (formaldehyde), Fix before staining (methanol), Fix after staining (methanol), Permeabilize after staining
Colors	Blue, Green, Orange, Red, Far-red, Near-infrared
Storage Conditions	See Product Information Sheet for details. Store at -10 to -35 °C. Protect fluorescent conjugates from light. After reconstitution store at 2-8 °C for up to 1 week, or at -10 to -35 °C for up to 12 months
Reconstitution	Dissolve 1 mg conjugate in 1 mL dH2O
Antibody/conjugate formulation	Lyophilized, 1 mg/mL in 1X PBS after reconstitution
Application Notes	Recommended staining concentration 1-5 ug/mL. See Product Information Sheet for detailed protocols
Shelf life	Guaranteed for at least 12 months from date of receipt when stored as recommended

Product	Conjugation	Ex/Em	Size	Catalog No.	Purchase
CF@350 WGA	CF@350	347/448 nm	1 mg 5 mg	29021-1 29021	Purchase 29021-1 Purchase 29021
CF@405S WGA	CF@405S	404/431 nm	1 mg 5 mg	29027-1 29027	Purchase 29027-1 Purchase 29027
CF@405M WGA	CF@405M	408/452 nm	1 mg 5 mg	29028-1 29028	Purchase 29028-1 Purchase 29028
CF@488A WGA	CF@488A	490/515 nm	1 mg 5 mg	29022-1 29022	Purchase 29022-1 Purchase 29022
CF@532 WGA	CF@532	527/558 nm	1 mg 5 mg	29064-1 29064	Purchase 29064-1 Purchase 29064
CF@555 WGA	CF@555	555/565 nm	5 mg	29076	Purchase 29076
CF@568 WGA	CF@568	562/583 nm	1 mg 5 mg	29076-1 29077	Purchase 29076-1 Purchase 29077
CF@594 WGA	CF@594	593/614 nm	1 mg 5 mg	29077-1 29023-1	Purchase 29077-1 Purchase 29023-1
CF@633 WGA	CF@633	630/650 nm	1 mg 5 mg	29023 29024-1	Purchase 29023 Purchase 29024-1
CF@640R WGA	CF@640R	642/662 nm	1 mg 5 mg	29024 29026-1	Purchase 29024 Purchase 29026-1
CF@680 WGA	CF@680	681/698 nm	1 mg 5 mg	29026 29029-1	Purchase 29026 Purchase 29029-1
CF@680R WGA	CF@680R	680/701 nm	1 mg 5 mg	29029 29025-1	Purchase 29029 Purchase 29025-1
CF@770 WGA	CF@770	770/797 nm	1 mg 5 mg	29025 29059-1	Purchase 29025 Purchase 29059-1
Biotin WGA	Biotin	N/A	1 mg 5 mg	29059 29095-1	Purchase 29059 Purchase 29095-1
HRP WGA	HRP	N/A	1 mg	29095 29073	Purchase 29095 Purchase 29073

This datasheet was generated on August 31, 2024 at 06:26:03 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/wheat-germ-agglutinin-wga-cf-dye-conjugates/>