

## Datura Stramonium Lectin (DSL) Conjugates



DSL binds to (beta-1,4) linked N-acetylglucosamine oligomers, preferring chitobiose or chitotriose over single N-acetylglucosamine residues. Biotium offers DSL conjugated to biotin and a selection of 6 bright and photostable CF® Dyes.

### Product Description

*Datura stramonium* lectin (DSL) binds to (beta-1,4) linked N-acetylglucosamine oligomers, preferring chitobiose or chitotriose over single N-acetylglucosamine residues. DSL also binds well to N-acetylglucosamine and oligomers containing repeating N-acetylglucosamine sequences. This lectin binds well in the acidic pH range, but its affinity decreases above pH 8.0. The lectin may be used to stain tissue sections. Biotium offers DSL conjugated to biotin and a selection of bright and photostable CF® Dyes.

- Binds to (beta-1,4) linked N-acetylglucosamine oligomers
- Suitable for immunofluorescence staining in tissue sections
- Choice of 7 CF® Dyes or biotin
- Supplied at 1 mg/mL in 10 mM HEPES pH 7.5, 0.15 M NaCl, 0.08% sodium azide, 0.1 mM CaCl<sub>2</sub>

**Note:** CF® Dye DSL conjugates are formulated with sodium azide and are not suitable for use *in vivo* or live cell culture

### Find the Right Stain for Your Application

DSL 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, [Ulex europaeus agglutinin I \(UEA I\)](#), [Lycopersicon esculentum \(Tomato\)](#), [Phaseolus vulgaris leucoagglutinin \(PHA-L\)](#), [Wheat Germ Agglutinin \(WGA\) conjugates](#), [Concanavalin A \(Con A\)](#) and [Arachis hypogaea \(PNA\)](#) 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 other fluorescent dyes. Learn more about [CF® Dyes](#).

### Datura Stramonium Lectin (DSL) Conjugates

Conjugation	Ex/Em	Size	Catalog No.	Dye Features
<a href="#">CF@488A</a>	490/516 nm	1 mL	<a href="#">29097</a>	<a href="#">CF@488A Features</a>
<a href="#">CF@568</a>	562/584 nm	1 mL	<a href="#">29098</a>	<a href="#">CF@568 Features</a>
<a href="#">CF@594</a>	593/615 nm	1 mL	<a href="#">29099</a>	<a href="#">CF@594 Features</a>
<a href="#">CF@640R</a>	642/663 nm	1 mL	<a href="#">29100</a>	<a href="#">CF@640R Features</a>
<a href="#">CF@680</a>	681/698 nm	1 mL	<a href="#">29101</a>	<a href="#">CF@680 Features</a>
<a href="#">CF@740</a>	742/767 nm	1 mL	<a href="#">29131</a>	<a href="#">CF@740 Features</a>
<a href="#">CF@790</a>	783/808 nm	1 mL	<a href="#">29146</a>	<a href="#">CF@790 Features</a>
<a href="#">Biotin</a>	N/A	1 mL	<a href="#">29096</a>	

### Full List of Lectin Conjugates

Product	Features
<a href="#">CF® Dye Concanavalin A (Con A)</a>	<ul style="list-style-type: none"> <li>• Cell surface stain for yeast, fungi, and mammalian cells</li> <li>• Selectively binds to α-mannopyranosyl and α-glucopyranosyl residues</li> <li>• Available with a wide selection of CF® Dyes</li> </ul>

### Product attributes

<b>Probe cellular localization</b>	Membrane/cell surface
<b>For live or fixed cells</b>	For fixed cells, For live/intact cells
<b>Cell permeability</b>	Membrane impermeant
<b>Fixation options</b>	Fix before staining (formaldehyde), Fix after staining (formaldehyde), Fix before staining (methanol), Fix after staining (methanol), Permeabilize after staining
<b>Colors</b>	Green, Red, Far-red, Near-infrared

Product	Features
<a href="#">CF® Dye Wheat Germ Agglutinin (WGA)</a>	<ul style="list-style-type: none"> <li>• Cell surface stain for mammalian cells and gram+ bacteria</li> <li>• Also stains yeast bud scars</li> <li>• Has high affinity for sialic acid and N-acetylglucosamine</li> <li>• Choose from a wide selection of CF® Dyes or HRP</li> </ul>
<a href="#">CF® Dye Peanut Lectin (PNA) from Arachis hypogaea</a>	<ul style="list-style-type: none"> <li>• Specific for terminal <math>\beta</math>-galactose and binds preferentially to galactosyl (<math>\beta</math>-1,3) N-acetylgalactosamine</li> <li>• Choice of 6 CF® dye colors</li> </ul>
<a href="#">CF® Dye Lycopersicon Esculentum (Tomato) Lectin (LEL, TL)</a>	<ul style="list-style-type: none"> <li>• Marker for blood vessels and microglial cells</li> <li>• Binds to [GlcNAc] 1,3-N-acetylglucosamine, glycophorin, and Tamm-Horsfall glycoprotein</li> <li>• Used to study tumor angiogenesis or tracing neovascular development in xenograft models</li> <li>• Choice of 7 CF® Dyes or biotin</li> </ul>
<a href="#">CF® Dye Ulex Europaeus Agglutinin I (UEA I)</a>	<ul style="list-style-type: none"> <li>• Marker for human endothelial cells and incompletely differentiated gastrin cells</li> <li>• Binds to glycoproteins and glycolipids containing <math>\alpha</math>-linked fucose residues</li> <li>• Choice of 7 CF® Dyes or biotin</li> </ul>
<a href="#">CF® Dye Phaseolus Vulgaris Leucoagglutinin (PHA-L)</a>	<ul style="list-style-type: none"> <li>• Used to stimulate lymphocyte and T cell proliferation</li> <li>• Choice of 7 CF® Dyes or biotin</li> </ul>
<a href="#">CF® Dye Datura Stramonium Lectin (DSL)</a>	<ul style="list-style-type: none"> <li>• Binds to (beta-1,4) linked N-acetylglucosamine oligomers</li> <li>• Choice of 7 CF® Dyes or biotin</li> </ul>
<a href="#">CF® Dye Sambucus Nigra Lectin (SNA, EBL)</a>	<ul style="list-style-type: none"> <li>• Binds to sialic acid attached to terminal galactose</li> <li>• Choice of 7 CF® Dyes or biotin</li> </ul>

CF is a registered trademark of Biotium, Inc.

This datasheet was generated on June 14, 2026 at 09:00:53 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/datura-stramonium-lectin-dsl/>