



# LECTIN KIT I, BIOTINYLATED

**SKU:** BK-1000



## DESCRIPTION

Our lectin screening kits are designed to provide the investigator with a panel of seven lectins or lectin conjugates. The lectins have been selected to offer a variety of sugar specificities and are of the same high quality as the reagents offered individually.

Each kit consists of 1 mg of the following lectins or lectin conjugates: Con A, DBA, PNA, RCA120, SBA, UEA I, WGA.

## SPECIFICATIONS

|                          |                                                                        |
|--------------------------|------------------------------------------------------------------------|
| <b>Unit Size</b>         | 1 Kit                                                                  |
| <b>Sugar Specificity</b> | Mannose, Galactose, Fucose, N-Acetylgalactosamine, N-Acetylglucosamine |

**For research use only. Not intended for therapeutic or diagnostic use in animals or humans.**



|                     |                                                                                                                      |
|---------------------|----------------------------------------------------------------------------------------------------------------------|
| <b>Applications</b> | Immunohistochemistry / Immunocytochemistry, Immunofluorescence, Blotting Applications, Elispot, ELISAs, Glycobiology |
| <b>Conjugate</b>    | Biotinylated                                                                                                         |

## KIT COMPONENTS

### Kit Contents:

This lectin screening kit is designed to provide the investigator with a panel of seven biotinylated lectin conjugates. The lectins have been selected to offer a variety of sugar specificities and are of the same high quality as the reagents offered individually.

Kit I consists of 1 mg of the following biotinylated lectins:

- Con A (Concanavalin A; 2 mg/ml)
- SBA (*Glycine max* (soybean) agglutinin; 2 mg/ml)
- WGA (*Triticum vulgaris* (wheat germ) agglutinin; 2 mg/ml)
- DBA (*Dolichos biflorus* agglutinin; 2 mg/ml)
- UEA I (*Ulex europaeus* agglutinin I; 2 mg/ml)
- RCA<sub>120</sub> (*Ricinus communis* agglutinin; 2 mg/ml)
- PNA (*Arachis hypogaea* (peanut) agglutinin; 2 mg/ml)

## TECHNICAL INFORMATION

These bioinylated conjugates are prepared from our affinity-purified lectins and are optimally labeled with biotin. Essentially free of inactive lectin conjugates and containing no free biotin, these biotinylated lectins provide ideal intermediates for examining glycoconjugates using the Biotin-Avidin/Streptavidin System. First a biotin-labeled lectin is added, followed by the VECTASTAIN ABC Reagent, Avidin D conjugate, or streptavidin derivative.

**For research use only. Not intended for therapeutic or diagnostic use in animals or humans.**



## Lectin Screening Kits

| Lectin                                      | Common Abbreviation | Primary Sugar Specificity        |
|---------------------------------------------|---------------------|----------------------------------|
| <b>Lectin Kit I</b>                         |                     |                                  |
| Concanavalin A                              | Con A               | Mannose                          |
| <i>Dolichos biflorus</i> agglutinin         | DBA                 | N-Acetylgalactosamine            |
| Peanut Agglutinin                           | PNA                 | Galactose                        |
| <i>Ricinus communis</i> agglutinin I        | RCA I               | Galactose, N-Acetylgalactosamine |
| Soybean agglutinin                          | SBA                 | N-Acetylgalactosamine            |
| <i>Ulex Europaeus</i> agglutinin I          | UEA I               | Fucose                           |
| Wheat Germ agglutinin                       | WGA                 | N-Acetylglucosamine              |
| <b>Lectin Kit II</b>                        |                     |                                  |
| <i>Griffonia simplicifolia</i> lectin I     | GSL I               | Galactose                        |
| <i>Len culinaris</i> lectin                 | LCA                 | Mannose                          |
| <i>Phaseolus vulgaris</i> Erythroagglutinin | PHA-E               | Complex structures               |
| <i>Phaseolus vulgaris</i> Leucoagglutinin   | PHA-L               | Complex structures               |
| <i>Pisum sativum</i> agglutinin             | PSA                 | Mannose                          |
| Wheat Germ agglutinin, succinylated         | Succinylated WGA    | N-Acetylglucosamine              |
| <i>Sophora japonica</i> agglutinin          | SJA                 | N-Acetylgalactosamine            |
| <b>Lectin Kit III</b>                       |                     |                                  |
| <i>Datura stramonium</i> lectin             | DSL                 | N-Acetylglucosamine              |
| <i>Erythrina cristagalli</i> lectin         | ECL                 | Galactose                        |
| <i>Griffonia simplicifolia</i> lectin II    | GSL II              | N-Acetylglucosamine              |
| Jacalin                                     | Jacalin             | Galactose                        |
| <i>Lycopersicon esculentum</i> lectin       | LEL                 | N-Acetylglucosamine              |
| <i>Solanum tuberosum</i> lectin             | STL                 | N-Acetylglucosamine              |
| <i>Vicia villosa</i> lectin                 | VVL                 | N-Acetylgalactosamine            |

## CITATIONS



Powered by Bioz © 2023 See more details on Bioz

## DOCUMENTS

- [Lectins in Histochemistry, ELISA, and Western Blot Applications](#)
- [Safety Data Sheet](#)
- [Download CoA](#)
- [Datasheet](#)

**For research use only. Not intended for therapeutic or diagnostic use in animals or humans.**



[www.vectorlabs.com](http://www.vectorlabs.com)

Email: [customerservice@vectorlabs.com](mailto:customerservice@vectorlabs.com)

Telephone: [\(650\) 697-3600](tel:(650)697-3600)

## GALLERY IMAGES



**For research use only. Not intended for therapeutic or diagnostic use in animals or humans.**

[Lectin Kit I, Biotinylated](https://vectorlabs.com/products/lectin-kit-i-biotinylated/)

<https://vectorlabs.com/products/lectin-kit-i-biotinylated/>