



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

Unit Size	1 Kit
Sugar Specificity	Mannose, Galactose, Fucose, N-Acetylgalactosamine, N-Acetylglucosamine

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



Applications	Immunohistochemistry / Immunocytochemistry, Immunofluorescence, Blotting Applications, Elispot, ELISAs, Glycobiology
Conjugate	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)
- RCA₁₂₀ (*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
Lectin Kit I		
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
Lectin Kit II		
<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
Lectin Kit III		
<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

Email: 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/>