



LECTIN KIT I, FLUORESCIN

SKU: FLK-2100



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 I (BK-1000, FLK-2100, RLK-2200) consists of 1 mg of the following lectins or lectin conjugates: Con A, DBA, PNA, RCA120, SBA, UEA I, WGA.

SPECIFICATIONS

Color of Fluorescence	Green
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	Immunofluorescence, Glycobiology
Conjugate	Fluorescein

KIT COMPONENTS

This lectin screening kit is designed to provide the investigator with a panel of seven fluorescein labeled 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 fluorescein labeled lectin conjugates:

- Con A (Concanavalin A; 2 mg/ml)
- SBA (*Glycine max* (soybean) agglutinin; 2 mg/ml)
- WGA (*Triticum vulgare* (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)

These lectins are supplied essentially free of unconjugated fluorochrome and inactive lectin.

Fluorescein labeled lectin has an excitation maximum at 495 nm and an emission maximum at 515 nm.

TECHNICAL INFORMATION

Our fluorescein labeled lectins are produced by using the highest quality fluorescein isothiocyanate, our affinity-purified lectins, and special conjugation procedures. The lectin conjugates have the appropriate number of fluorochromes bound which provide the maximum fluorescence and optimum staining characteristics for each particular lectin. These lectins are supplied essentially free of unconjugated fluorochrome and inactive lectin. Fluorescein labeled lectin has an excitation maximum at 495 nm and an emission maximum at 515 nm.

Accompanying each fluorescent lectin kit is an analysis data sheet summarizing the results of our quality control tests and providing pertinent information on the product. All of these reagents are supplied as solutions preserved with sodium azide.

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, Fluorescein](https://vectorlabs.com/products/lectin-kit-i-fluorescein)

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