



LECTIN KIT II, BIOTINYLATED

SKU: BK-2100



DESCRIPTION

Our lectin screening kits are designed to provide the investigator with a panel of six 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 II (LK-3000, BK-2100, FLK-3100, RLK-3200) consists of 1 mg of the following lectins or lectin conjugates: GSL I, LCA, PHA-E, PHA-L, PSA, Succinylated WGA.

SPECIFICATIONS

Unit Size	1 Kit
Applications	Immunohistochemistry / Immunocytochemistry, Immunofluorescence, Blotting Applications, Elispot, ELISAs, Glycobiology
Conjugate	Biotinylated

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



KIT COMPONENTS

This lectin screening kit is designed to provide the investigator with a panel of six 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 II consists of 1 mg of the following biotinylated lectins:

- GSL I (*Griffonia (Bandeiraea) simplicifolia* lectin I; 2 mg/ml)
- PSA (*Pisum sativum* agglutinin; 2 mg/ml)
- LCA (*Lens culinaris* agglutinin; 2 mg/ml)
- PHA-E (*Phaseolus vulgaris* Erythroagglutinin; 2 mg/ml)
- PHA-L (*Phaseolus vulgaris* Leucoagglutinin; 2 mg/ml)
- Succinylated WGA (Wheat germ agglutinin, succinylated; 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.

DOCUMENTS

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

GALLERY IMAGES

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)



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

[Lectin Kit II, Biotinylated](https://vectorlabs.com/products/lectin-kit-2-biotinylated/)

<https://vectorlabs.com/products/lectin-kit-2-biotinylated/>