



# Lectin Kit I, Biotinylated

## BK-1000

Product Images



## Short 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.

## Additional Information

---

Unit Size	1 kit
Applications	Immunohistochemistry / Immunocytochemistry, Immunofluorescence, Blotting Applications, Elispot, ELISAs, Glycobiology
Conjugate	Biotinylated
Sugar Specificity	Mannose, Galactose, Fucose, N-Acetylgalactosamine, N-Acetylglucosamine

