



# Lectin Kit III, Biotinylated

## BK-3000

Product Images



## Additional Information

---

Unit Size	1 kit
Applications	Immunohistochemistry / Immunocytochemistry, Immunofluorescence, Blotting Applications, Elispot, ELISAs, Glycobiology
Recommended Usage	Reconstitute lyophilized biotinylated lectins by adding 0.5 ml water. The resulting solution will be: 10 mM HEPES, 0.15 M NaCl, 0.1 mM CaCl <sub>2</sub> , 0.08% sodium azide, pH 7.5.
Conjugate	Biotinylated
Sugar Specificity	Galactose, N-Acetylgalactosamine, N-Acetylglucosamine

