



# Streptavidin/Biotin Blocking Kit SP-2002

Product Images



## Short Description

---

Streptavidin/Biotin Blocking Kit blocks all endogenous biotin, biotin receptors, and streptavidin binding sites present in tissues. This kit is designed for use with streptavidin detection systems such as the VECTASTAIN® Universal Quick kits if streptavidin or biotinylated reagents bind non-specifically to tissues or proteins. This blocking kit consists of 18 ml Streptavidin and 18 ml of biotin in convenient dropper bottles.

## Additional Information

---

Unit Size	1 kit
Applications	Immunohistochemistry / Immunocytochemistry, Immunofluorescence, In situ hybridization, Blotting Applications, Elispot
Blocking Action	Endogenous Biotin or Avidin/Streptavidin Binding Proteins
Format	Ready-to-Use

