



# Avidin/Biotin Blocking Kit

## SP-2001

Product Images



## Short Description

---

Avidin/Biotin Blocking Kit blocks all endogenous biotin, biotin receptors, and avidin binding sites present in tissues. This kit is designed for use with avidin detection systems such as the VECTASTAIN® ABC Kits if avidin or biotinylated reagents bind non-specifically to tissues or proteins.

This blocking kit consists of 18 ml of Avidin D 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

