



# **AVIDIN/BIOTIN BLOCKING KIT**

**SKU:** SP-2001



---

## **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.

## **SPECIFICATIONS**

**Blocking Action** Endogenous Biotin or Avidin/Streptavidin Binding Proteins

**Format** Ready-to-Use

**Unit Size** 1 Kit

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



## Applications

Immunohistochemistry / Immunocytochemistry, Immunofluorescence,  
In situ hybridization, Blotting Applications, Elispot

# TECHNICAL INFORMATION

Product review on [Biocompare.com](http://Biocompare.com).

## CITATIONS



Powered by Bioz © 2023 See more details on Bioz

## DOCUMENTS

- [User Guide](#)
- [Safety Data Sheet](#)
- [Download CoA](#)
- [Datasheet](#)

## GALLERY IMAGES



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