



STREPTAVIDIN/BIOTIN BLOCKING KIT

SKU: SP-2002



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.

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

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.