



HORSERADISH PEROXIDASE AVIDIN D (AV-HRP), CONCENTRATE, FOR ELISAS AND BLOTS

SKU: A-2014-5



DESCRIPTION

Vector Laboratories' enzyme-conjugated avidin and streptavidin are produced with the highest specific activity enzymes in optimal ratios. Specific covalent linkages are chosen to provide stable, highly active conjugates.

Horseradish peroxidase labeled Avidin D (Av-HRP), for ELISAs and Blots, is optimized for use in solid-phase assays such as ELISAs and blotting applications.

SPECIFICATIONS

Detection Enzyme	Peroxidase
Format	Concentrate
Formulation	0.05 M NaHCO ₃ , pH 8.3, preservative

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



Unit Size	5 mg
Storage Instructions	2-8 °C DO NOT FREEZE
Usage Summary	For diluting this product, we recommend a buffered saline solution, approximately pH 7.5. The recommended concentration range for use is 2-5 µg/ml.
Applications	Blotting Applications, Elispot, ELISAs
Concentration	5 mg active conjugate/ml
Conjugate	Peroxidase

TECHNICAL INFORMATION

For HRP Avidin conjugate that is optimized for tissue/cell staining systems please see the following product:

[Avidin D, Horseradish Peroxidase \(Av-HRP\), Concentrate, for IHC](#)

DOCUMENTS

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

GALLERY IMAGES



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