



Horseradish Peroxidase Avidin D (Av-HRP), Concentrate, for ELISAs and Blots

A-2014-5

[Product Images](#)



Short 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.

Additional Information

Unit Size	5 mg
Applications	Blotting Applications, Elispot, ELISAs
Recommended Usage	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.
Solution	0.05 M NaHCO ₃ , pH 8.3, preservative
Recommended Storage	2-8 °C DO NOT FREEZE
Concentration	5 mg active conjugate/ml
Detection Enzyme(s)	Peroxidase
Conjugate	Peroxidase
Format	Concentrate

