



STREPTAVIDIN, PEROXIDASE, ELISA GRADE (CONCENTRATE, FOR ELISAS AND BLOTS)

SKU: SA-5014-1



DESCRIPTION

Features:

- Avidin and streptavidin reagents are powerful tools to detect or purify biotinylated proteins, nucleic acids, and other macromolecules.
- Optimized for use in solid phase assays such as ELISAs and blotting applications
- Horseradish Peroxidase Streptavidin (HRP Streptavidin) is produced by our own coupling procedure which preserves the high specific activity of the peroxidase
- For horseradish peroxidase streptavidin (Sav-HRP) that is optimized for use in tissue/cell staining applications, see the following product:

Streptavidin, Horseradish Peroxidase (Sav-HRP), Concentrate, for IHC/ICC

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



SPECIFICATIONS

| | |
|-----------------------------|--|
| Detection Enzyme | Peroxidase |
| Format | Concentrate |
| Formulation | 0.05 M HEPES, 0.1M Na2S04pH 6.8, preservative |
| Unit Size | 1 mg |
| Storage Instructions | 2-8 °C, DO NOT FREEZE |
| Usage Summary | Do not use sodium azide with this product since it is an inhibitor of horseradish peroxidase. Recommended concentration range for use is 5-10 µg/ml. Note: This product is different from Cat. No. SA-5004 which has been optimized for use in immunohistochemical applications. |
| Applications | Blotting Applications, Elispot, ELISAs |
| Concentration | 1.0 mg/ml active conjugate |
| Conjugate | Peroxidase |

TECHNICAL INFORMATION

Avidin and streptavidin reagents are powerful tools to detect or purify biotinylated proteins, nucleic acids, and other macromolecules.

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.

This Horseradish peroxidase streptavidin preparation (HRP-Sav; SA-5014) is optimized for use in solid-phase assays such as ELISAs and blotting applications.

For horseradish peroxidase streptavidin (Sav-HRP) that is optimized for use in tissue/cell staining applications, see the following product:

[Streptavidin, Horseradish Peroxidase \(Sav-HRP\), Concentrate, for IHC/ICC](#)

CITATIONS

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



Powered by Bioz © 2023 See more details on Bioz

DOCUMENTS

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

GALLERY IMAGES



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