



Streptavidin, Peroxidase, ELISA Grade (Concentrate, for ELISAs and Blots)

SA-5014-1

[Product Images](#)



Short 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](#)

Additional Information

Unit Size	1 mg
Applications	Blotting Applications, Elispot, ELISAs
Recommended Usage	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.
Solution	0.05 M HEPES, 0.1M Na ₂ SO ₄ pH 6.8, preservative
Recommended Storage	2-8 °C, DO NOT FREEZE
Concentration	1.0 mg/ml active conjugate
Detection Enzyme(s)	Peroxidase
Conjugate	Peroxidase
Format	Concentrate

