
Streptavidin: HRP Conjugate

(Cat. No. R-1098)

Description

Streptavidin: horseradish peroxidase conjugates show negligible non-specific binding to non-biotinylated macromolecules and therefore give very low backgrounds.

Purification

Purified Streptavidin conjugated to Horseradish Peroxidase (HRP) - liquid

Concentration

1 mg/ml

Formulation

Phosphate buffered saline, 40% glycerol

Specificity

Biotinylated macromolecules

Storage

Store at -20°C. This product should be stored undiluted.

Handling Recommendations

Should this product contain a precipitate we recommend microcentrifugation before use. For maximum recovery of the products, centrifuge the vial prior to opening the cap.

Applications & Suggested Dilutions

ELISA:1:1000 – 1:5000; Immunohistology - Frozen:1:100-1:500; Immunohistology - Paraffin:1:200; Where this product has not been tested for use in a particular technique this does not necessarily exclude its use in such procedures. Suggested working dilutions are given as a guide only. It is recommended that the user titrates the protein for use in their own system using appropriate negative/positive controls.

Ordering Information

Products	Size	Cat. No.
Streptavidin: HRP Conjugate	0.5 ml	R-1098-1

This product is for research purposes only. Not intended for use in diagnostic procedures.