

Product Information

SigmaScreen™ Streptavidin coated plates, 96 Well Plates

Catalog Numbers:

M5432	Clear
M5557	White opaque
M4058	Black opaque
M3433	Clear strips

Storage Temperature 2-8 °C

TECHNICAL BULLETIN

Product Description

Streptavidin multiwell plates are coated with streptavidin, a 60 kDa protein isolated from *Streptomyces avidinii*. The purified protein is bound to the wells of polystyrene plates via a proprietary coating technology. This coating technology ensures:

- high binding capacity of biotin
- high coating homogeneity
- low leaching of streptavidin (<6 ng per well)
- high resistance to commonly used detergents

In addition, streptavidin-coated multiwell plates are pre-blocked for immediate use.

Binding assays of biotinylated single and double stranded DNA, peptides, proteins, and small organic molecules can be performed on streptavidin-coated multiwell plates.

Reaction Volume

Streptavidin is coated at a reaction volume of 100 µl/well. The wells are blocked at 200 µl/well.

Binding Capacity

In saturation and competitive binding assays performed on this product, binding of ≥5 pmoles of d-biotin per well is observed. The binding capacity for larger molecules labeled with biotin may be less than that of d-biotin due to steric hindrance associated with the specific molecule.

Precautions and Disclaimer

This product is for R&D use only, not for drug, diagnostic procedures, household, or other uses. Please consult the Material Safety Data Sheet for information regarding hazards and safe handling practices.

Storage/Stability

For optimal performance, the unopened product should be stored in a dry place at 2–8 °C. Under these storage conditions, the product is stable for two years. For short-term storage of less than 3 months, the product may be stored at room temperature. Once opened, it is suggested that the product be used immediately.

Not recommended for assays at >60 °C.

Procedures

Plate Viability Assay

To validate the viability of the streptavidin surface follow the procedure below:

1. Dissolve 1 mg of biotinylated horseradish peroxidase (P9568) in 1 ml of Phosphate buffered saline (PBS) containing 0.05% TWEEN® 20 (P3563). Dilute the 1 mg/ml stock 1:50,000–1:100,000 in PBS containing 0.05% TWEEN 20 and add 100 µl per well. As a negative control, add 100 µl of the same dilution of streptavidin-peroxidase (S5512) to a separate set of wells.
2. Incubate the wells for 30 minutes at room temperature.
3. Wash the wells three times, 300 µl per well, with PBS containing 0.05% TWEEN 20.
4. After discarding the final wash, add 100 µl per well of TMB substrate (T8665).
5. Incubate the wells for 15 minutes before reading the absorbance in a spectrophotometer. If desired, the reaction may be stopped with the addition of 0.5 M H₂SO₄ (50 µl per well). An absorbance of at least 1.5 will be observed at 655 nm for a non-stopped reaction or at 450 nm for a stopped reaction.

Peptide and Protein Binding

1. Prepare a solution of the biotinylated protein or peptide in either PBS or Tris buffered saline (TBS), pH 7.4. A starting concentration of 1–10 µg/ml should be used if the optimal concentration is not known.
2. Add up to 100 µl of the solution per well and allow the samples to incubate for 1–2 hours at a temperature within the range of 18–30 °C. Include blank and control wells as appropriate.
3. Wash the wells three times, 300 µl per well, with PBS or TBS containing 0.05% TWEEN 20 (P3563 or T9039).
4. Incubate the wells with 100 µl of an appropriately diluted primary antibody in PBS or TBS containing 0.05% TWEEN 20 for 30 minutes to 1 hour.
5. Wash the wells three times, 300 µl per well, with PBS or TBS containing 0.05% TWEEN 20.
6. Incubate the wells with 100 µl of an appropriately diluted enzyme-labeled secondary antibody in PBS or TBS containing 0.05% TWEEN 20 for 30 minutes to 1 hour.
7. Wash the wells three times, 300 µl per well, with PBS or TBS containing 0.05% TWEEN 20.
8. After addition of an appropriate substrate, the wells are ready for detection by various modes (colorimetry, chemiluminescence or fluorescence).

PCR Products

1. Prepare the biotinylated PCR product for addition onto streptavidin-coated multiwell plates by diluting the sample 1:10-1:50 in PBS containing 0.05% TWEEN 20.
2. Apply 100 µl per well of the diluted sample and allow the sample to incubate for 30–60 minutes at 25–37 °C.
3. Bound PCR products are denatured by adding 0.5 M NaOH, 100 µl per well. Incubate for 5–10 minutes.
4. Wash the wells three times, 300 µl per well, with PBS containing 0.05% TWEEN 20 to remove the nonbiotinylated, complimentary strand of the PCR product.
5. Add 200 µl per well of a hapten-labeled oligonucleotide that is complimentary to the biotinylated strand. Use 0.05–0.5 pmole of labeled oligonucleotide per well. Hybridize in the presence of 5× SSC buffer with 0.3% TWEEN 20 and 1% BSA. Allow the hybridization to proceed for 30–60 minutes at 37–50 °C.
6. Wash the wells three times, 300 µl per well, with PBS containing 0.05% TWEEN 20.
7. Add 100 µl per well of an appropriately diluted detection conjugate in PBS containing 0.05% TWEEN 20. Incubate for 30–60 minutes at room temperature.
8. Wash the wells five to six times, 300 µl per well, with PBS containing 0.05% TWEEN 20.
9. Detect the bound PCR products using an appropriate substrate (see below).

Results

Optimization of ELISA Results

There are four major areas where detection can be optimized: nonspecific binding, wash conditions, antibody affinity, and conjugate concentration.

1. Nonspecific Binding:

Factors that contribute to nonspecific binding are ionic interactions, hydrophobic interactions, and cross-reactivity. To reduce nonspecific binding, changes in conjugate concentrations and wash buffers can be made. Users are encouraged to modify buffers with components in the concentration ranges indicated below.

Detergents	0.05–0.1% TWEEN 20 (P9416) 0.02–0.1% CHAPS (C5070)
Salt	0.5–1.0 M NaCl (S3014) or Na ₂ HPO ₄ (S3264)
Protein blockers	0.1–1% BSA (A9647), Casein (C3400), or Gelatin (G7765)
Non protein blockers	1% PEG 20 (P2263) or Polyvinylpyrrolidone (P5288)

2. Wash Conditions:

To limit reversible nonspecific binding interactions, at least three wash steps are recommended.

3. Antibodies and Conjugates:

For optimal signal performance, the user is encouraged to use high-affinity antibodies and conjugates. Commercially obtained antibodies and conjugates should be used at the concentrations suggested by the supplier.

References

1. Wilchek, M., and Bayer, E.A., (eds.), Avidin-Biotin Technology. *Meth. Enzymol.*, **184** (1990).
2. Crowther, J.R. (ed.), ELISA: Theory and Practice. *Methods in Molecular Biology*, Vol **42** (Humana Press, Totowa, NJ, 1995), Product code Z36,415-0.
3. Dekonenko, A., *et al.*, A colorimetric PCR-enzyme immunoassay to identify Hantaviruses. *Clin and Diagnostic Virol.*, **8**, 113-121 (1997).
4. Morozov, I., *et al.*, Mapping of functional domains in p47^{phox} involved in the activation of NADPH oxidase by "peptide walking". *J. Biol. Chem.*, **273**, 15435-15444 (1998).
5. Hemminki, A., *et al.*, Fine tuning of an anti-testosterone antibody binding site by stepwise optimisation of the CDRs. *Immunotechnol.*, **4**, 59-69 (1998).
6. Rong, R., *et al.*, Two-site immunofluorometric assay of intact salmon calcitonin with improved sensitivity. *Clin. Chem.*, **43**, 71-75 (1997).
7. Vesanen, M., *et al.*, Detection of Herpes simplex virus DNA in cerebrospinal fluid samples using the polymerase chain reaction and microplate hybridization. *J. Virol. Meth.*, **59**, 1-11 (1996).
8. Benotmane, A.M., *et al.*, Nonisotopic quantitative analysis of protein-DNA interactions at equilibrium. *Anal. Biochem.*, **250**, 181-185 (1997).
9. Verhaegen, M., and Christopoulos, T.K., Quantitative polymerase chain reaction based on a dual-analyte chemiluminescence hybridization assay for target DNA and internal standard. *Anal. Chem.*, **70**, 4120-4125 (1998)

CDP-*Star* is a registered trademark of Tropix, Inc. Bedford, MA, USA, and covered under US patent 5,326,882.

SigmaScreen is a trademark of Sigma-Aldrich Biotechnology LP and Sigma-Aldrich Co.

TWEEN is a registered trademark of Croda International PLC

RG,PHC 07/10-1

Sigma brand products are sold through Sigma-Aldrich, Inc.

Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications.

Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply.

Please see reverse side of the invoice or packing slip.