



SIGMA[®]

**Multiwell Plate, 96 Wells
Streptavidin Coated**

Product Numbers: Z382469, Z382477, Z382485

MULTIWELL PLATE, 96 WELLS STREPTAVIDIN COATED

Technical Bulletin

INTENDED USE

For Research Use Only
Not for use in diagnostic procedures.

PRODUCT DESCRIPTION

Streptavidin multiwell plates are coated with streptavidin, a 60 KD protein isolated from *Streptomyces avidinii*. The purified protein is bound to the wells of polystyrene microtiter 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.

High capacity 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 μ l/well. The wells are blocked at 200 μ l/well.

BINDING CAPACITY

In saturation and competitive binding assays performed on this product, binding of ≥ 25 pmoles of d-biotin per well is observed. The binding capacity of larger molecules labeled with biotin will possibly be less than that of biotin via steric factors associated with the specific molecule.

STORAGE AND HANDLING CONDITIONS:

For optimal performance, the unopened product should be stored in a dry place at 2–30°C. Under these storage conditions, the product is stable for two years. Once opened, it is suggested that the product be used immediately.

Not recommended for assays at $> 60^{\circ}\text{C}$.

SAMPLE PROTOCOLS FOR THE USE OF STREPTAVIDIN COATED MULTIWELL PLATES:

PLATE VIABILITY ASSAY

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

1. Using catalog item P 9568 (Biotinylated HRP), dissolve 1 mg in 1 ml of PBS + 0.05% Tween-20 (P 3563).
Dilute the 1 mg/ml stock 1:50,000–1:100,000 in PBS + 0.05% Tween-20 and add 100 μ l per well. As a negative control, add 100 μ l of a similar dilution of streptavidin-peroxidase (S 5512) to a separate set of wells.
2. At room temperature, incubate the wells for 30 minutes.

3. Wash the wells three times, 300 μ l per well, with PBS + 0.05% Tween-20.
4. After discarding the final wash, add 100 μ l per well of TMB substrate (T 8665).
5. Incubate the wells for 15 minutes before reading in a spectrophotometer. If desired, the reaction may be stopped with the addition of 0.5M H_2SO_4 (50 μ l per well). An absorbance of at least 1.5 will be observed at 655 nm for a non-stopped reaction or 450 nm for a stopped reaction.

PEPTIDE AND PROTEIN BINDING

1. Prepare a solution of the biotinylated protein or peptide in 0.05M bicarbonate buffer pH 9.6, PBS pH 7.4 or TBS pH 7.5. A starting concentration of 1–10 μ g/ml should be used if the optimal concentration is not known.

2. Add up to 100 μ l of the solution per well and allow the samples to incubate for 1–2 hours at a temperature range of 18–30°C.
3. Wash the wells three times, 300 μ l per well, with PBS or TBS + 0.05% Tween-20.
4. Incubate the wells with 100 μ l of an appropriately diluted primary antibody in PBS or TBS + 0.05% Tween-20 for 30 minutes to 1 hour.
5. Wash the wells three times, 300 μ l per well, with PBS or TBS + 0.05% Tween-20.
6. Incubate the wells with 100 μ l of an appropriately diluted enzyme labeled secondary antibody in PBS or TBS + 0.05% Tween-20 for 30 minutes to 1 hour.
7. Wash the wells three times, 300 μ l per well, with PBS or TBS + 0.05% Tween-20.
8. After addition of an appropriate substrate, the wells are ready for detection in various modes (colorimetry, chemiluminescence or fluorescence).

PCR PRODUCTS

1. Prepare the biotinylated PCR product for addition onto streptavidin coated multiwell plates by diluting the reaction 1:10–1:50 in PBS + 0.05% Tween-20.
2. Apply 100 μ l per well of the diluted product and allow the sample to incubate for 30–60 minutes at 25–37°C.
3. Bound PCR products are denatured by adding 100 μ l per well of 0.5M NaOH. Incubate for 5–10 minutes.
4. Wash the wells three times, 300 μ l per well, with PBS + 0.05% Tween-20 to remove the nonbiotinylated, complimentary strand of the PCR product.
5. Add 200 μ l per well of a hapten labeled oligonucleotide which is complimentary to the biotinylated strand. Use 0.05–0.5 pmole of labeled oligonucleotide per well. Hybridize in

the presence of 5X SSC, 0.3% Tween-20, 1% BSA. Allow the hybridization to proceed for 30–60 minutes at 37–50°C.

6. Wash the wells three times, 300 μ l per well, with PBS + 0.05% Tween-20.
7. Add 100 μ l per well of an appropriately diluted detection conjugate in PBS + 0.05% Tween-20. Incubate for 30–60 minutes at RT.
8. Wash the wells five to six times, 300 μ l per well, with PBS + 0.05% Tween-20.
9. Proceed to the use of the specific substrate for detection.

DETECTION

There exists a number of alternatives for the detection of molecules labeled with common haptens. Below is a table of indirect detection systems for two haptens, namely biotin and fluorescein.

	Common Colorimetric Substrates	Common Chemiluminescent Substrates and Activators
Peroxidase	TMB, ABTS	Luminol/ Iodophenol/peroxide
Alkaline phosphatase	pNPP	dioxetanes
Acridinium ester		NaOH/H ₂ O ₂
Beta-galactosidase		dioxetanes

OPTIMIZATION OF ELISA RESULTS

There are four major areas where detection results can be optimized: nonspecific binding, antibody affinity, wash conditions, 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 ranges of the various conditions stated below.

Detergents	0.05–0.1% Tween-20 0.002–0.05% Tween-65
Salt	0.5–1.0 M NaCl or Na ₂ HPO ₄
Protein blockers	0.1–1% BSA or casein
Non protein blockers	1% PEG 20 or Polyvinylpyrrolidone

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.

RELATED PRODUCTS

- T 8665 3,3',5,5'-Tetramethylbenzidine (TMB) Liquid Substrate System
- A 9941 2,2'-Azino-bis(3-ethylbenzthiazoline-6-sulfonic acid) tablets
- P 4922 Phosphate-citrate buffer tablets with Sodium Perborate capsules
- N 7653 p-Nitrophenyl Phosphate (pNPP) Liquid Substrate System

- P 3813 Phosphate Buffered Saline, pH 7.4
- P 3688 Phosphate Buffered Saline with BSA, pH 7.4
- P 3563 Phosphate Buffered Saline with Tween-20, pH 7.4
- T 6664 Tris Buffered Saline, pH 8.0
- T 9039 Tris Buffered Saline with Tween-20, pH 8.0
- P 1318 Phosphatase, Alkaline-biotinamidocaproyl labeled
- P 9568 Peroxidase-biotinamidocaproyl labeled
- S 5512 Streptavidin-peroxidase labeled
- S 6639 20X SSC Buffer concentrate
- A 8327 30% BSA solution
- P 9416 Polyoxyethylenesorbitan monolaurate (Tween-20)
- BK-200 Cleavable Biotinylation Kit
- BK-101 Immunoprobe Biotinylation Kit
- B-TAG Biotin Tag Micro Biotinylation Kit
- OLB-16-DDU BioProbe 3'Oligonucleotide Labeling Kit with Bio-16-ddUTP

GENERAL REFERENCES

1. Wilchek M, Bayer EA: eds, Meth Enzymol, Avidin-Biotin Technology, 184, 1990
2. Crowther JR: ed, "ELISA: Theory and Practice" Methods in Molecular Biology, Vol 42 (Humana Press, Totowa, NJ, 1995), Catalog # Z36,415-0
3. Dekonenko A, Ibrahim MS, Schmaljohn CS: A Colorimetric PCR-Enzyme Immunoassay to Identify Hantaviruses. Clin and Diagnostic Virology, 8:113-121, 1997
4. Fujita S: *et al*, Microtitration Plate Enzyme Immunoassay To Detect PCR-Amplified DNA from Candida Species in Blood. J Clin Microbiol, 33:962-967, 1995
5. Guenther PC, Hart CE: Quantitative, Competitive PCR Assay for HIV-1 Using a Microplate-Based Detection System. Biotechniques, 24:810-816, 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 AM: *et al*, Nonisotopic Quantitative Analysis of Protein-DNA Interactions at Equilibrium. Anal Biochem, 250:181-185, 1997
9. Verhaegen M, Christopoulos TK: Quantitative Polymerase Chain Reaction Based on a

Dual-Analyte Chemiluminescence Hybridization Assay for Target DNA and Internal Standard. Anal Chem, 70:4120-4125, 1998

Sigma-Aldrich, Inc. warrants that its product 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 invoice or packing slip.

Previous Revision: April 1999
Revised: September 1999

SIGMA CHEMICAL COMPANY

P.O. BOX 14508, ST. LOUIS, MO 63178 USA

In the USA/Canada call: **800-325-3010** for orders and **800-325-8070** for customer service
Outside the USA/Canada call: 314-286-7750 for orders and 314-289-8470 for customer service