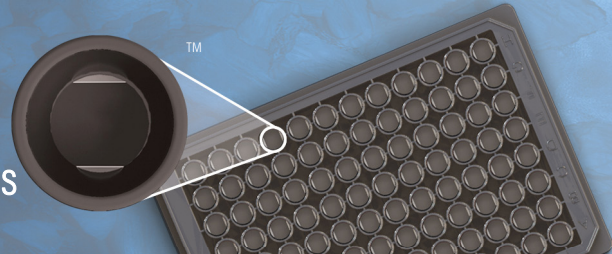


New Product

MULTI-ARRAY Streptavidin and High Bind Avidin Gold Plates



Catalog Numbers

MULTI-ARRAY 96-well Streptavidin Gold Plates for SECTOR [®] Imager	
Kit size	
1 plate	L15SA-1
5 plates	L15SA-2
30 plates	L15SA-5

MULTI-ARRAY 96-well High Bind Avidin Gold Plates for SECTOR Imager	
Kit size	
1 plate	L15AB-1
5 plates	L15AB-2
30 plates	L15AB-5

MULTI-ARRAY 96-well Streptavidin Gold Plates for SECTOR PR [®]	
Kit size	
1 plate	L13SA-1
5 plates	L13SA-2
30 plates	L13SA-5

Visit www.mesoscale.com for a complete listing of available materials and reagents.

Ordering information

MSD Customer Service
 Phone: 1-301-947-2085
 Fax: 1-301-990-2776
 Email: CustomerService@mesoscale.com

Company Address

MESO SCALE DISCOVERY
 A division of
 Meso Scale Diagnostics, LLC.
 9238 Gaither Road
 Gaithersburg, MD 20877 U.S.A.

MESO SCALE DISCOVERY[®] launches a new gold standard in streptavidin and high bind avidin-coated plates ensuring unsurpassed performance, consistency and convenience for immunogenicity testing, pharmacokinetic/pharmacodynamic studies, and assay development.

Key Advantages

- Improved Plate Performance:** Stringent specifications set for consistency, uniformity, and binding capacity
- Guaranteed Lot-to-lot Consistency:** Every lot meets consistent release criteria as specified on the full Certificate of Analysis
- Increased Shelf Life:** 30-month shelf life of Streptavidin Gold Plates assures performance in longitudinal studies
- Large Plate Lots:** Standard Streptavidin Gold Plate lot sizes of 3000, larger lots available on request
- Assured Availability:** Orders shipped within 48 hours of order acceptance

The Certificate of Analysis included with each shipment confirms the tighter specifications.

Sample Certificate of Analysis

Product Description: MULTI-ARRAY[®] 96-well Streptavidin Gold Plate
Catalog #: L15SA-1 / 1 plate; L15SA-2 / 5 plates; L15SA-5 / 30 plates; L15SA-6 / 120 plates; L15SA-7 / 510 plates
Lot Number: _____
Expiration date: _____
Statement: This product was manufactured and tested according to approved MSD[®] Standard Operating Procedures.

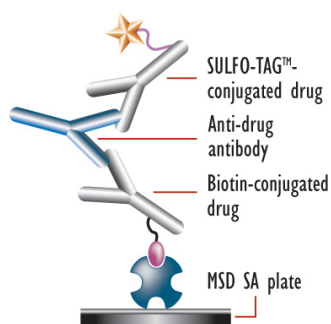
Streptavidin Whole Plate Uniformity QC Specifications		
Metric	Specification	Pass/Fail?
Mean Intraplate CV	≤ 6%	Pass
Intraplate CVs	≤ 8%	Pass
Number of plates with intraplate CV >12%	0 plates	Pass
Interplate CV	≤ 8%	Pass
Plates where signal > 20% from plate mean occurs in same well on multiple plates	0 plates	Pass
Wells with signal > 50% from plate mean	0 plates	Pass
Median signal for concentric rings, max to min range	≤ 10% for all plates	Pass
Median signal for columns, max to min range	≤ 10%	Pass
Median signal for rows, max to min range	≤ 10%	Pass
Number of plates sampled (N)	Dependent on Lot Size	Pass

Binding Capacity QC Specifications			
Concentration of SULFO-TAG [™] Labeled Biotinylated IgG (BTI)	Mean ECL specification and tolerance range	Intraplate %CV or SD specification	Pass/Fail?
6 nM (0.3 pmole)	82,211 counts ± 15%	10%	Pass
4 nM (0.2 pmole)	55,593 counts ± 15%	10%	Pass
2 nM (0.1 pmole)	26,330 counts ± 15%	10%	Pass
0 nM	≤ 100 counts	15 count SD	Pass

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

Bridging Immunogenicity Assay on the MSD Platform

MSD electrochemiluminescence has been the choice for the development of immunogenicity assays since 2003. MSD assays offer multiple advantages over common ELISAs.



Desirable Assay Requirements	ELISA	MSD
Free Drug Tolerance	Poor	Excellent
Detection of Low Affinity Antibodies	No	Yes
Few Wash Steps	3-4	1
High Throughput	Good	High
Direct Conjugation of Stable Label	Yes	Yes
High Level of Sensitivity	100s ng/mL	10s ng/mL
Wide Dynamic Range	1-2 logs	3-4 logs
Minimal Sample Volume	25-100 µL	5-25 µL
High Binding Capacity		10x more

Example Assay Protocol: The protocol is simple, reproducible and scalable.

1. Combine biotin-drug, SULFO-TAG drug and sample in a polypropylene plate. Incubate for 1 hour to overnight.
2. Transfer solution to pre-blocked MULTI-ARRAY Streptavidin Gold Plate. Incubate for 1 hour.
3. Wash. Add Read Buffer. Read plate.

Additional MSD Advantages

- Flexible assay formats enable immunogenicity testing of many drug types, including antibodies, proteins, and peptides
- Rapid assay development yields efficient and cost effective results
- On-site technical support by MSD's field-based Application Scientists is available for seminars, training and assistance with assay development
- Leading contract research organizations (CROs) are using MSD's technology to run validated assays

MSD is committed to meeting the ever-increasing demands of the immunogenicity community. Call us at 240-631-2522 or visit our website at www.mesoscale.com and put us to work for you.