

B-068 MULTI-ARRAY[®] Assay to Discriminate Recent from Long-Standing HIV Infection

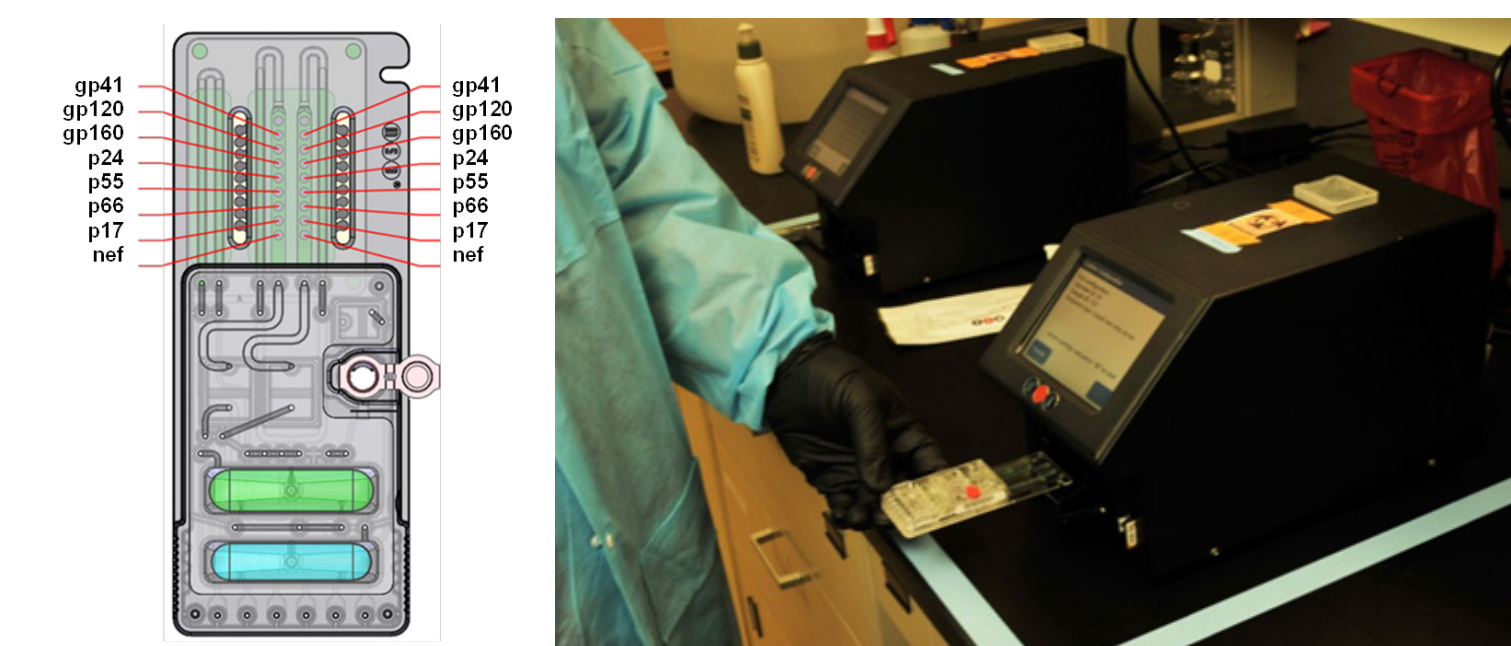
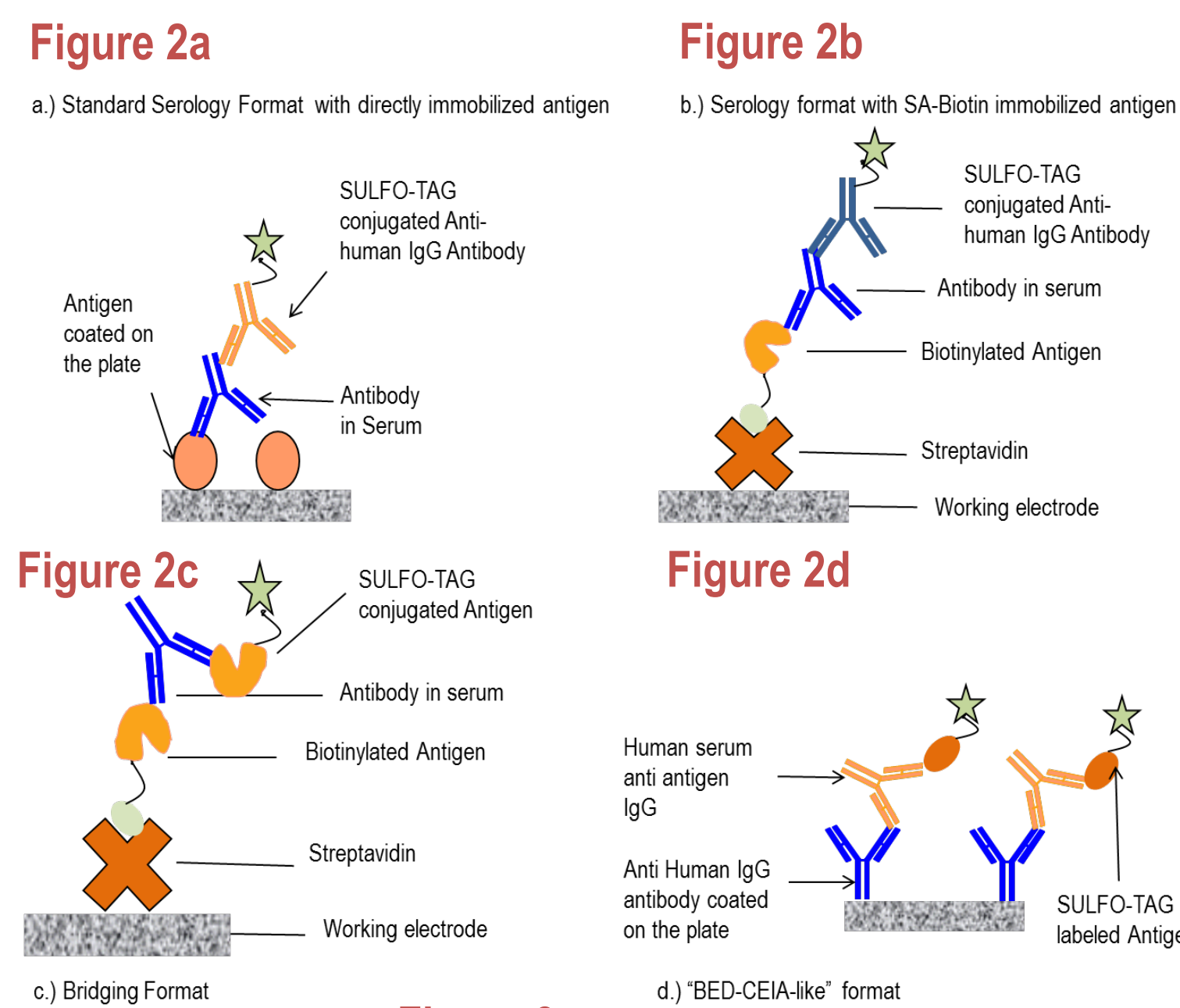
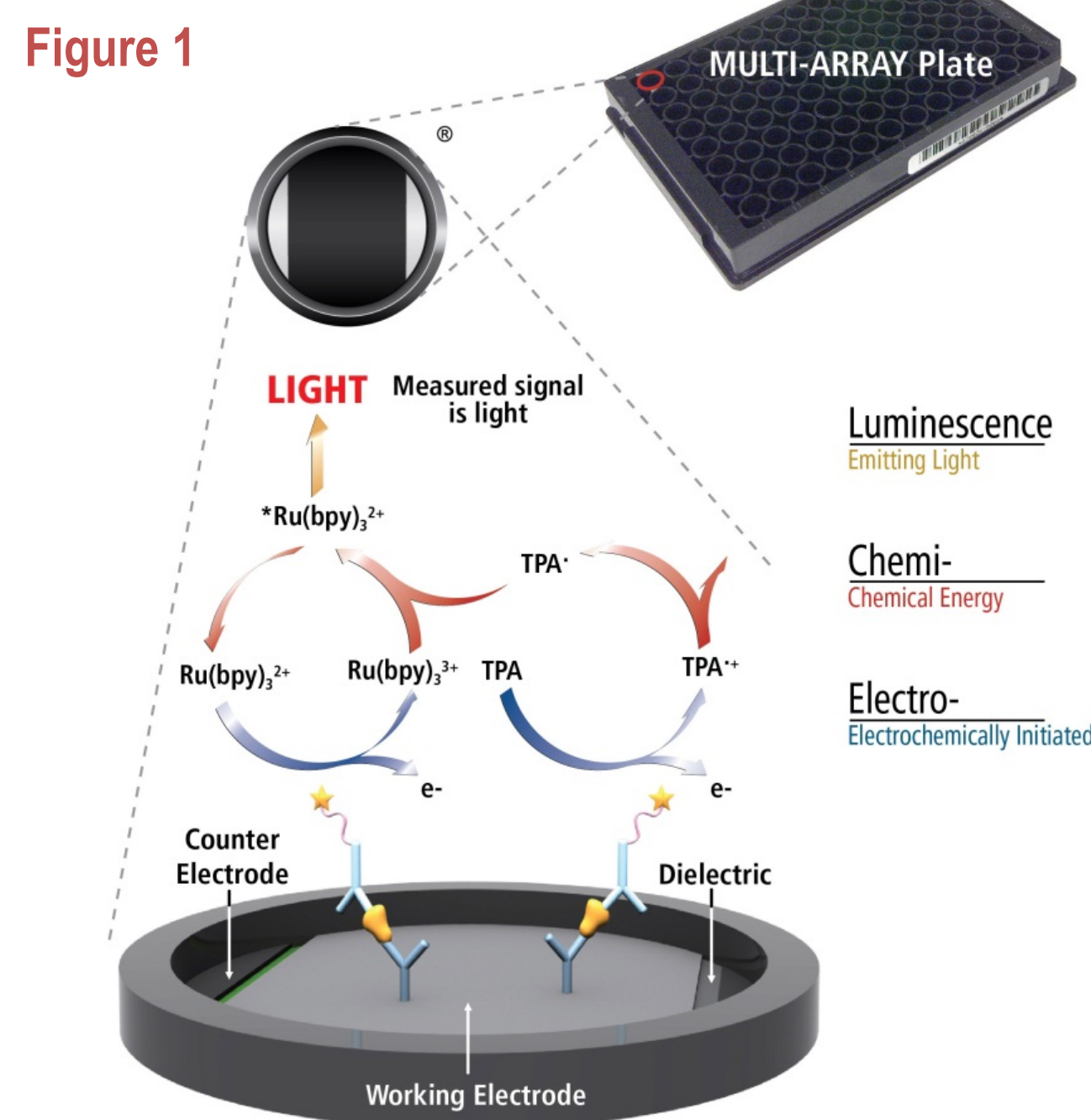
Martin Stengelin, Daisy Roy, Mikayla Higgins, Iva Maxwell, Zara Karuman, Eli N. Glezer, and Jacob N. Wohlstadter
Meso Scale Diagnostics, LLC. Rockville, Maryland, USA

1 Abstract

In order to accurately assess and compare different prevention strategies, the rate at which new HIV infections are acquired in a population needs to be measured accurately. A simple laboratory test that indicates whether an HIV infection was acquired in the recent past (generally 4-12 months) would be very useful to estimate HIV incidence. We demonstrated feasibility of several assay formats to separate recent from longstanding HIV infection. Using MULTI-ARRAY technology, we measured antibodies against the HIV proteins gp41, gp120, gp160, p17, p24, p55, p66, tat, viv, and nef in a multiplexed format using a very small sample volume (25 μ L of a 1,000-fold diluted serum or plasma sample). We used the well-characterized "HIV Incidence/Prevalence Performance Panel" from SeraCare (part # PRB601), which contains plasma samples from 15 HIV positive donors that have been characterized either as "incident" (recent infection) or "prevalent" (longstanding infection) based on consensus results from nine tests. Our MULTI-ARRAY serology format for antibodies against gp120 and gp160 showed ~10-fold separation between the median signals for incident and prevalent samples (and another ~10-fold separation from apparently healthy controls). All samples in each of the three groups were completely separated from the other two groups. We also developed avidity assay formats for antibodies against gp41, gp120, gp160, and p66 that could accurately separate samples from patients with incident versus prevalent HIV infection. The assays were developed in a 96-well high-throughput assay format for the MESO[®] SECTOR S 600 Imager and the MESO QuickPlex[®] SQ 120. We demonstrated feasibility for transfer of the assay format to a point-of-care (POC) platform. The POC assay is fully automated and simultaneously measures concentrations of antibodies against eight HIV proteins. Time to result is 25 minutes, and CVs are approximately 13%. The magnitude of the antibody response against gp120 and against gp160 accurately separates patients with incident HIV infection from patients with prevalent HIV infection, equivalent to the plate-based results. In conclusion, we demonstrated feasibility for development of high-throughput and point-of-care assays to discriminate recent from longstanding HIV infection.

2 Methods

MSD's electrochemiluminescence detection technology uses SULFO-TAG[™] labels that emit light upon electrochemical stimulation initiated at the electrode surfaces of MULTI-ARRAY and MULTI-SPOT[®] microplates.



POC Platform

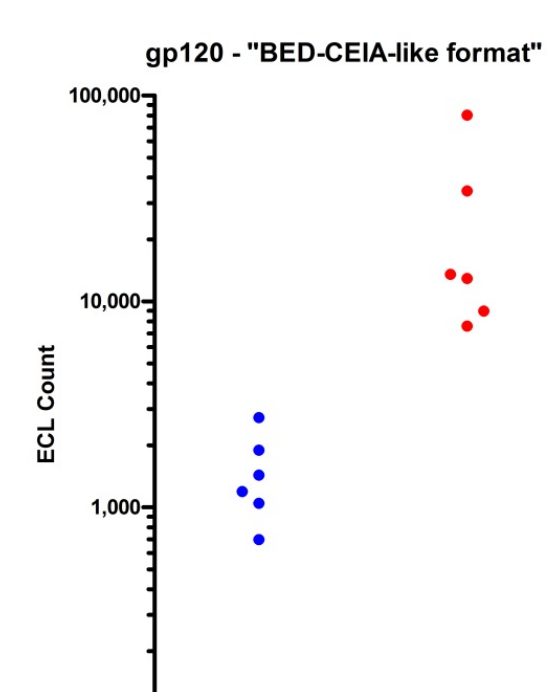
Figure 3 shows the MSD POC single-use cartridge and the cartridge reader. One sample is processed per cartridge; however, the cartridge has two independent detection channels that can be run with the same sample but using different reagent formulations, providing some flexibility for supporting assays with different optimal processing conditions. Two patterned arrays of antigens (or antibodies) are contained in the cartridge, immobilized on screen-printed carbon ink electrodes in the two independent detection channels.

3 MULTI-ARRAY Bridging Assay Format and "BED-CEIA-Like" Format

Feasibility of the bridging format in a single-plex mode was demonstrated using biotinylated antigen and streptavidin coated plates. The data on the right for 125-fold diluted plasma samples from the SeraCare Incident/Prevalent panel show a clear difference between disease states for all four tested HIV antigens.

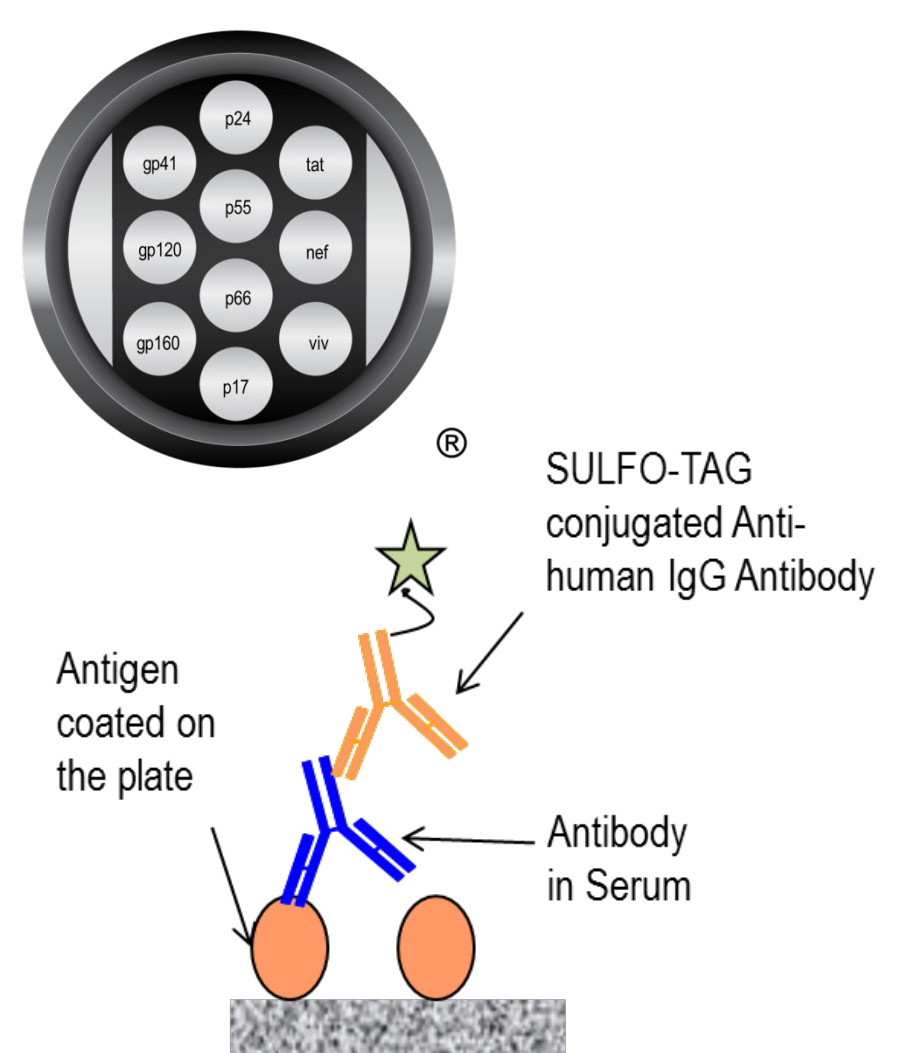
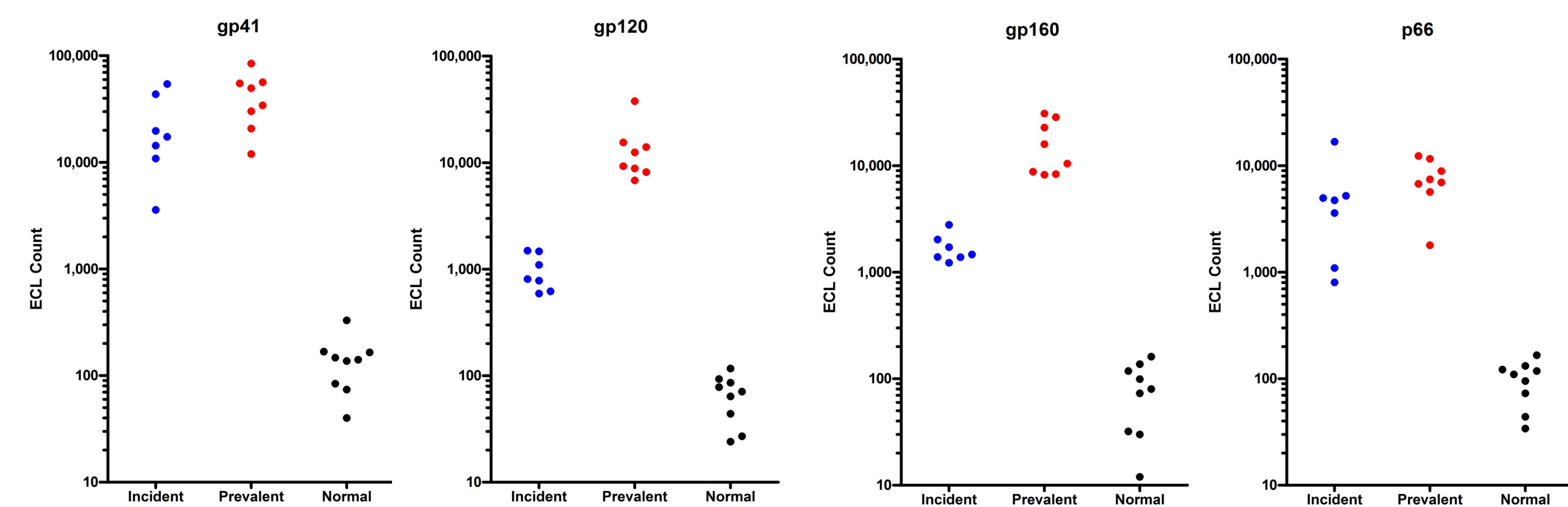
	Bridging Format: ECL Counts				
	gp41	gp120	gp160	p24	
Incident	PRB601-01	16,880	7,089	6,512	7,842
	PRB601-02	13,774	3,270	4,255	9,139
	PRB601-05	5,180	9,962	1,849	7,146
	PRB601-07	41,886	4,606	9,084	2,004
Prevalent	PRB601-03	48,630	76,019	29,294	216,253
	PRB601-04	41,296	144,517	21,137	108,286
	PRB601-06	49,855	93,393	25,083	1,742
	PRB601-08	54,104	125,977	10,533	192,696
Normal Samples	Donor-1	886	537	404	451
	Donor-2	947	543	358	465
	Donor-3	792	480	393	492
	Donor-4	483	551	379	442

Plates were coated with anti-human IgG, and after sample addition SULFO-TAG labeled HIV antigen (e.g. gp120) was added. This format allows discriminating incident from prevalent HIV infection.



4 Plate-Based MULTI-ARRAY Direct Serology Assay

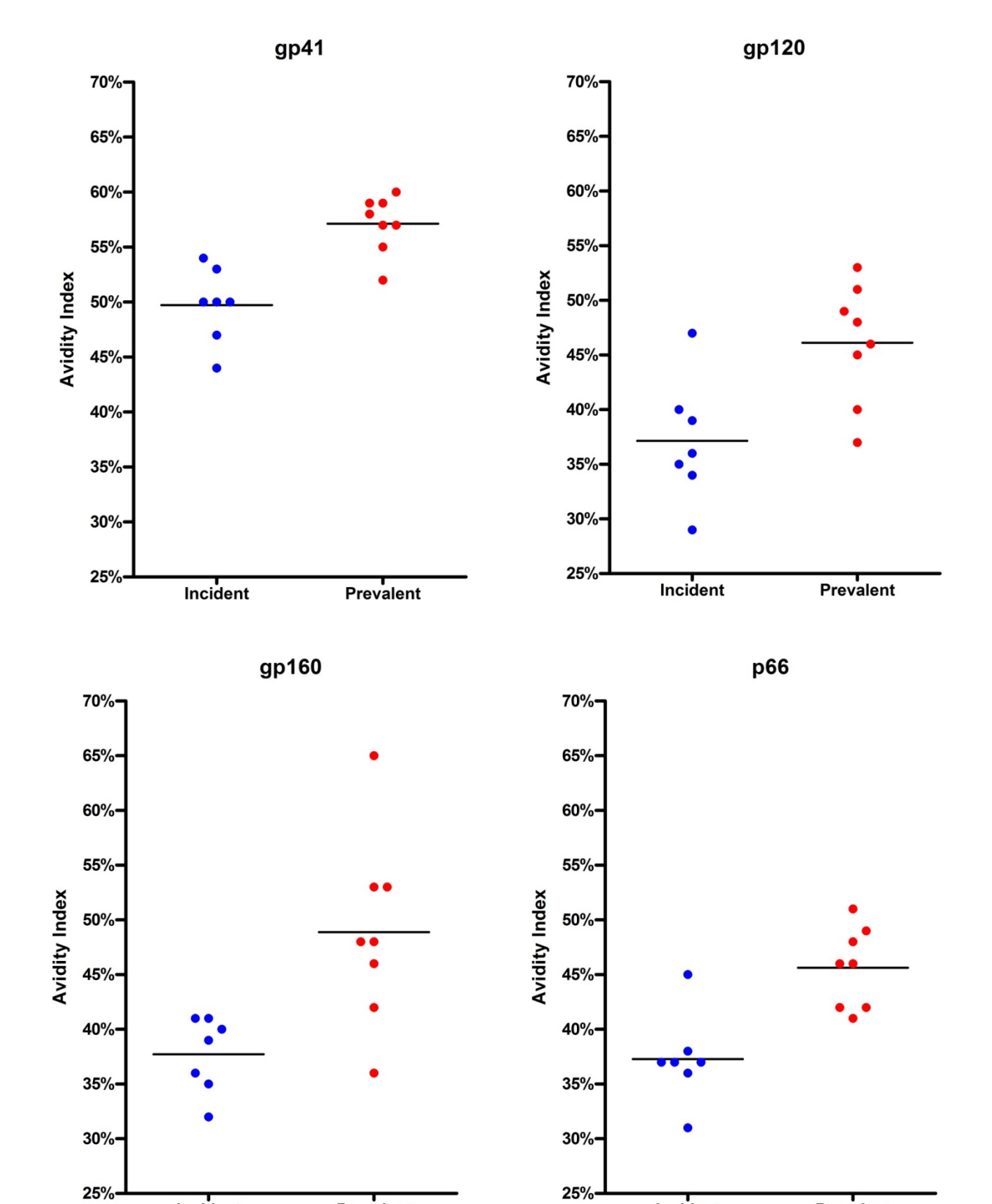
MULTI-ARRAY Serology Assay: ECL Counts for 1,000-fold diluted Plasma Sample											
Sample ID	Incident / Prevalent	gp41	gp120	gp 160	p24	p55	p66	p17	tat	nef	viv
PRB601-01	Incident	17,336	1,470	2,782	10,387	144	4,966	246	212	1,109	465
PRB601-02	Incident	19,701	590	1,225	11,153	126	5,207	209	181	900	554
PRB601-05	Incident	14,336	1,489	1,468	8,401	129	1,092	120	122	681	257
PRB601-07	Incident	43,540	620	1,717	2,921	178	801	160	210	841	349
PRB601-09	Incident	54,248	779	2,024	10,619	132	3,594	175	138	1,285	286
PRB601-12	Incident	3,582	1,096	1,387	3,556	144	4,725	180	198	1,133	748
PRB601-14	Incident	10,847	807	1,381	2,661	135	16,753	127	125	458	491
PRB601-03	Prevalent	30,146	9,263	8,204	16,020	268	1,792	177	187	1,450	533
PRB601-04	Prevalent	20,723	13,985	8,335	4,346	159	6,752	189	137	149	1,228
PRB601-06	Prevalent	49,533	8,152	10,460	3,939	166	7,421	123	192	424	312
PRB601-08	Prevalent	54,983	37,798	28,488	72,773	1,282	8,901	3,751	417	524	501
PRB601-10	Prevalent	56,419	12,504	22,759	165	156	6,962	2,160	166	523	416
PRB601-11	Prevalent	34,240	6,815	15,908	348	118	12,329	136	220	4,805	428
PRB601-13	Prevalent	84,232	15,484	30,780	5,291	117	5,659	172	157	490	257
PRB601-15	Prevalent	11,972	8,838	8,752	26,909	154	11,597	165	105	777	279



5 Plate-Based MULTI-ARRAY Avidity Assay

MULTI-ARRAY Serology Assay: Avidity Index											
Sample ID	Incident / Prevalent	gp41	gp120	gp 160	p24	p55	p66	p17	tat	nef	viv
Incident PRB601-01	Incident	50%	39%	41%	43%	45%	36%	51%	39%	40%	49%
Incident PRB601-02	Incident	54%	47%	41%	56%	24%	38%	15%	26%	43%	62%
Incident PRB601-05	Incident	50%	29%	32%	38%	61%	31%	64%	70%	42%	61%
Incident PRB601-07	Incident	47%	34%	36%	31%	42%	37%	36%	22%	42%	34%
Incident PRB601-09	Incident	53%	36%	35%	31%	32%	37%	28%	49%	45%	36%
Incident PRB601-12	Incident	50%	40%	40%	50%	33%	37%	28%	21%	44%	33%
Incident PRB601-14	Incident	44%	35%	39%	27%	45%	45%	2%	34%	40%	47%
Prevalent PRB601-03	Prevalent	60%	49%	48%	56%	65%	46%	39%	64%	58%	43%
Prevalent PRB601-04	Prevalent	55%	51%	53%	53%	45%	51%	39%	52%	43%	48%
Prevalent PRB601-06	Prevalent	59%	40%	42%	30%	31%	42%	53%	36%	35%	50%
Prevalent PRB601-08	Prevalent	59%	53%	65%	53%	35%	49%	37%	36%	15%	34%
Prevalent PRB601-10	Prevalent	52%	48%	48%	10%	37%	46%	34%	17%	32%	16%
Prevalent PRB601-11	Prevalent	57%	45%	46%	41%	33%	48%	31%	15%	50%	20%
Prevalent PRB601-13	Prevalent	58%	46%	53%	39%	30%	41%	51%	8%	37%	42%
Prevalent PRB601-15	Prevalent	57%	37%	36%	60%	88%	42%	83%	99%	51%	62%

A MULTI-ARRAY serology assay was performed as in Section 4, but for each sample, an additional condition was tested: 0.5 M of GuHCl was added to the read buffer, followed by a 5-minute incubation. The ratio of the signals in the presence and absence of GuHCl was defined as the avidity index. The table above shows the avidity index for all ten HIV proteins for the SeraCare HIV Incidence/Prevalence Performance Panel. As the graphs on the right demonstrate, there was a significant difference in the avidity index for incident versus prevalent samples for gp41, gp120, gp160, and p66.

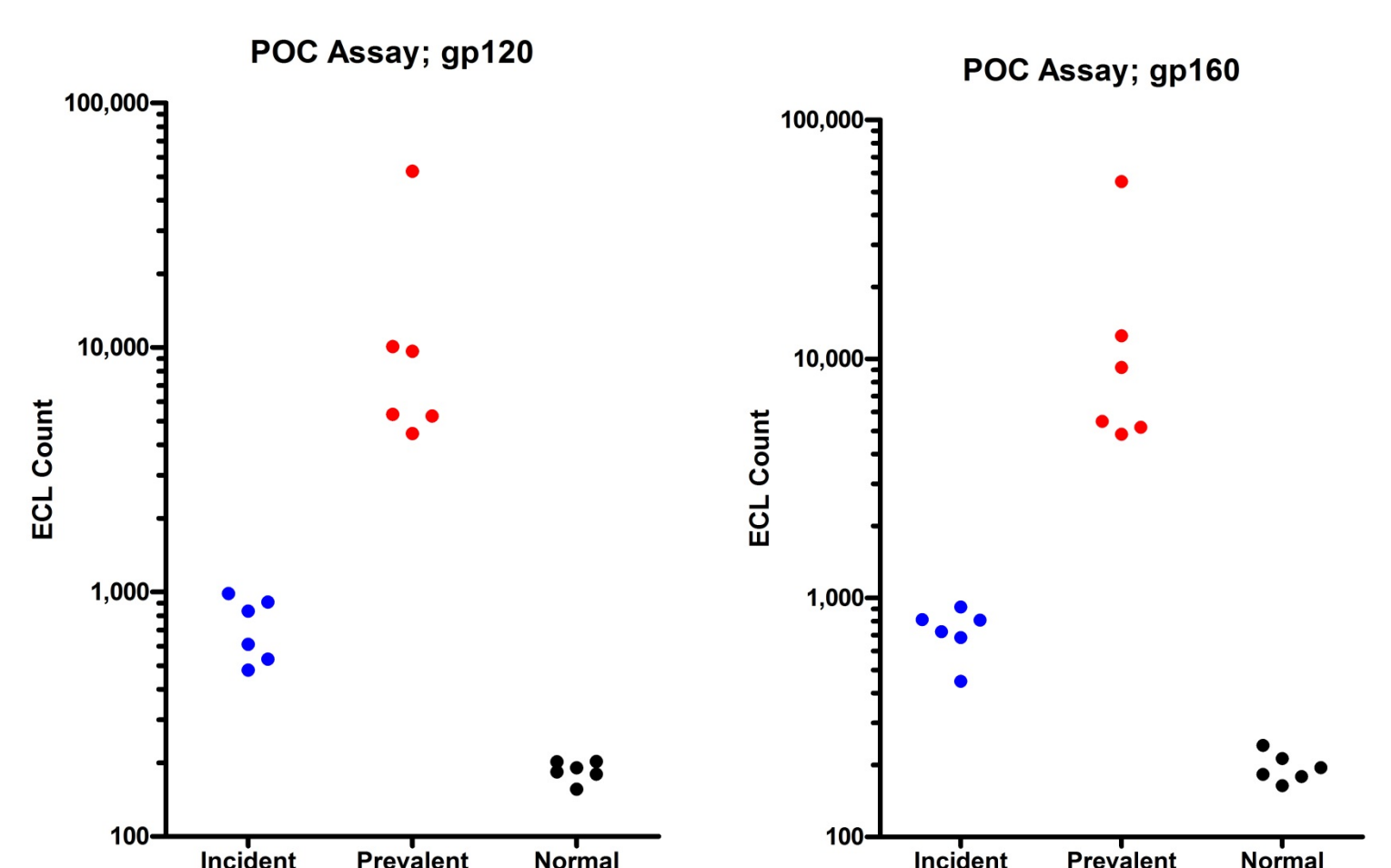


6 Feasibility of Transfer to POC Platform

The MSD POC cartridge includes two independent detection channels, each with eight spots. We immobilized recombinant gp41, gp120, gp160, p24, p55, p66, p17, and nef in both detection areas. SULFO-TAG labeled anti-human IgG detection antibody and buffer were dried in a dedicated area of the cartridge which can be hydrated and mixed by moving the sample back and forth. After a predetermined incubation time, the fluid is then moved to the capture spots for a second incubation followed by a wash and read.

The data on the right are from a 2-step assay format, where diluted sample was allowed to bind to the spots, followed by a wash and incubation with detection antibody. Time to result is 25 minutes, and CVs are approximately 13%.

In this sample set, the POC test could accurately differentiate all incident versus long-standing HIV infections, and could also resolve both groups from non-infected individuals.



7 Conclusion

We demonstrated feasibility for development of high-throughput and point-of-care assays to discriminate recent from longstanding HIV infection. Based on our results with the SeraCare HIV Incidence/Prevalence Performance Panel, the most promising markers to discriminate incident from prevalent HIV infection are the magnitudes of the antibody responses to gp120 and to gp160. Also promising are avidity to gp41, gp120, gp160, and p66.

As a next step, these results need to be confirmed with a larger set of clinical samples.

Acknowledgement:

The following reagents were obtained through the NIH AIDS Reagent Program, Division of AIDS, NIAID, NIH: HIV-1 nef Protein (Cat# 11478), HIV-1 viv Protein (Cat# 11050). This project has been funded in whole or in part with Federal funds from the National Institute of Health, Department of Human Health and Services, under Contract No. HHSN261201000018C.



Meso Scale Discovery[®]
A division of Meso Scale Diagnostics, LLC.
www.mesoscale.com[®]

DOWNLOAD POSTER



MESO SCALE DISCOVERY, MESO SCALE DIAGNOSTICS, MSD, MSD GOLD, DISCOVERY WORKBENCH, MULTI-ARRAY, MULTI-SPOT, QUICKPLEX, SECTOR, SECTOR PR, SECTOR HTS, SULFO-TAG, U-PLEX, S-PLEX, V-PLEX, STREPTAVIDIN GOLD, MESO, www.mesoscale.com, SMALL SPOT (design), 96 WELL 1, 4, 7 & 10-SPOT (design), 384 WELL 1 & 4-SPOT (design), MSD (design), U-PLEX (design), S-PLEX (design), V-PLEX (design), and SPOT THE DIFFERENCE are trademarks and/or service marks of Meso Scale Diagnostics, LLC. ©2015 Meso Scale Diagnostics, LLC. All rights reserved.