Serological Assays for Early Detection of HPV-associated Oropharyngeal Cancer

Mingyue Wang¹, John C. Smith¹, Jermaine Brown¹, Leonid Dzantiev¹, Yeming Wang¹, Anu Mathew¹, Martin Stengelin¹, George Sigal¹, Padhma Yuvaraj², Karen S. Anderson², and Jacob Wohlstadter¹ ¹Meso Scale Diagnostics, LLC. (MSD), Rockville, MD; ²Center for Personalized Diagnostics, Biodesign Institute, Arizona State University, Tempe, Arizona

1 Abstract

Research Use Only (RUO) sample-sparing MULTI-ARRAY serology immunoassays have been developed for human antibodies to several HPV-16 early antigens. These assays are intended to support research of oropharyngeal cancer (OPC) linked to the human papillomavirus (HPV). The HPV virus can infect the mouth and throat and cause cancers of the oropharynx, i.e. the back of the throat, including the base of the tongue and tonsils. HPV-related cancer incidence is increasing dramatically among men and is unlikely to decrease in the foreseeable future, despite vaccination efforts. As a result, this may become one of the most common cancers in middle-aged men in the United States by 2045. The HPV-16 genome consists of six early genes (E1, E2, E4, E5, E6, and E7) and two late genes (L1 and L2) that constitute the viral capsid. Antibodies against the late HPV-16 antigens are common in individuals with HPV infections and are observed in approximately 15% of the general population. Antibodies to early HPV-16 antigens are less common (about 1% of general population) but are commonly present in those with OPC. These antibodies are present many years before cancer develops; thus, they are ideal biomarkers for early detection of OPC. Immunoassays for antibodies against HPV-16 early antigens E1, E2, E6, and E7 were developed in a high-throughput multiplexed format using electrochemiluminescence (ECL) detection, and MULTI-ARRAY 10- spot 96-well plates. Each well has an array presenting recombinant antigens manufactured by MSD. The assay uses 25 uL of 2,500x diluted serum or plasma. The assay format is simple: diluted sample is incubated in the well with the antigen array followed by a wash and detection of bound anti-HPV antibodies using an anti-human IgG antibody labeled with the SULFO-TAG ECL label. Assay performance was evaluated with approximately 200 commercially sourced samples from apparently healthy individuals and 14 samples from individuals known to be positive for at least one HPV-16 early antigen using an established reference method (Programmable Protein (RAPID) ELISA, Anderson 2015). There was excellent agreement between the two methods. This high-throughput MULTI-ARRAY assay may be useful in research applications requiring screening of a large number of samples for HPV-16 antibodies.

2 Methods

MSD[®] 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.





MULTI-ARRAY Electrochemiluminescence Technology

- Minimal non-specific background and strong responses to analyte yield high signal-to-background ratios.
- The stimulation mechanism (electricity) is decoupled from the response (light signal), minimizing matrix interference. • Only labels bound near the electrode surface are excited, enabling non-
- washed assays. • Labels are stable, non-radioactive, and directly conjugated to biological
- molecules. • Emission at ~620 nm eliminates problems with color quenching. • Multiple rounds of label excitation and emission enhance light levels and
- improve sensitivity. Carbon electrode surface has 10X greater binding capacity than polystyrene wells.
- Surface coatings can be customized.

3 HPV-16 E-Protein Serology Assay

A multiplexed serology panel was developed to detect antibodies against HPV-16 E-proteins associated with oropharyngeal cancer. Plates are provided with antigens on spots in the wells of a 96-well plate. Antibodies in the sample bind to the antigens on the spots and anti-human IgG antibody conjugated with MSD SULFO-TAG is used for detection. Commercially-sourced plasma or serum samples were tested at 2,500-fold dilution on the serology panel.

	Anti-HPV E1 IgG	Anti-HPV E2 lgG	Anti-HPV E6 IgG	Anti-HPV E7 IgG
Positivity Rate in HPV-16 (+) Control Samples (N=14)	79%	93%	79%	93%
Positivity Rate in Breast Cancer Samples (N=40)	23%	23%	5%	18%
Positivity Rate in Gastric Cancer Samples (N=12)	25%	8%	5%	0%
Positivity Rate in Ovarian Cancer Samples (N=17)	0%	0%	6%	0%
Positivity Rate in Lung Cancer Samples (N=62)	7%	2%	15%	5%
Positivity Rate in Apparently Healthy (n=199)	5%	5%	5%	5%

Table 1. Sensitivity of seropositivity for cancer-associated HPV-16 E-Proteins at a 95% specificity.

Assay Protocol

- 1. Add diluent (25 μ L/per well) to the plates coated with antigens.
- 2. Add sample (25 µL/per well of 2,500-fold diluted serum or plasma).
- 3. Incubate 2 hours at room temperature (RT).
- 4. Wash and add detection antibody solution (25 µL per well). Incubate 1 hour at RT.
- 5. Wash and add read buffer (150 µL per well). Analyze with MSD instrument.

Individual Serum and Plasma Sample Testing

More than 300 commercially-sourced serum and plasma samples from apparently healthy individuals and from individuals with lung, breast, gastric or ovarian cancer, and 14 samples from individuals known to be positive for serological response to at least one HPV-16 early antigen were tested for HPV-16 E1, E2,E6, E7, HPV-16 L1 and HPV-18 L1 immunoglobulin G (IgG) antibodies. For oropharyngeal cancer-associated E proteins, preliminary cutoff values for positive serology were defined as the 95th percentile of signals from approximately 200 serum and plasma samples from apparently healthy individuals. Larger studies will be required to set a more clinically appropriate cutoff (e.g. 98-99th percentiles).

Age, gender and HPV vaccine status are not known for the samples tested in this study. We did not show a cutoff for L1 serology. A reasonable estimate for expected seropositivity for HPV-16 L1 in the general population is ~15%.



Figure 1. Absolute ECL signals for assays detecting antibodies against HPV-16 E proteins, HPV-16 L1, and HPV-18 L1 in commercially-sourced samples. For E-proteins, dashed black lines show the 95th percentile of signals from approximately 200 serum and plasma samples from apparently healthy individuals.





6 Results: Samples Show Positive Reactivity Against HPV-16 Early Antigen

Sample ID	Anti-HPV E1 IgG	Anti-HPV E2 lgG	Anti-HPV E6 IgG	Anti-HPV E7 IgG	Number of Pos Reactivities
HPV-16 Ab (+) Control 01	Pos	Pos	Neg	Pos	3
HPV-16 Ab (+) Control 02	Pos	Pos	Pos	Pos	4
HPV-16 Ab (+) Control 03	Pos	Pos	Pos	Pos	4
HPV-16 Ab (+) Control 04	Neg	Pos	Pos	Pos	3
HPV-16 Ab (+) Control 05	Pos	Pos	Neg	Pos	3
HPV-16 Ab (+) Control 06	Pos	Pos	Neg	Neg	2
HPV-16 AD (+) Control U/	Pos	POS	Pos	Pos	4
HPV-16 AD (+) Control 00	Pos	Pos	POS	POS	4 2
HPV-16 Ab (+) Control 10	Pos	Pos	Pos	<u> </u>	Δ
HPV-16 Ab (+) Control 11	Pos	Pos	Pos	Pos	4
HPV-16 Ab (+) Control 12	Pos	Pos	Pos	Pos	4
HPV-16 Ab (+) Control 13	Neg	Pos	Pos	Pos	3
HPV-16 Ab (+) Control 14	Pos	Pos	Pos	Pos	4
Breast Cancer 02	Neg	Neg	Neg	Pos	1
Breast Cancer 03	Pos	Pos	Neg	Pos	3
Breast Cancer 05	Neg	Pos	Neg	Neg	1
Breast Cancer 08	Pos	Pos	Neg	Neg	2
Breast Cancer 09	Pos	Neg	Neg	Neg	1
Breast Cancer 10	Pos	Pos	Pos	Pos	4
Breast Cancer 11	Pos	Pos	Neg	Pos	3
Breast Cancer 12	Pos	Pos	Neg	Pos	3
Breast Cancer 13	Pos	Neg	Neg	Neg	1
Breast Cancer 14	POS	POS	Neg	POS	3
Broast Cancer 10	Reg	Pos	Neg	Neg	
Gastric Cancer 04	Pos	Neg	Neg	Neg	1
Gastric Cancer 07	Pos	Pos	Neg	Neg	2
Gastric Cancer 12	Pos	Neg	Neg	Neg	1
Ovarian Cancer 12	Neg	Neg	Pos	Neg	1
Lung Cancer 11	Pos	Neg	Pos	Neg	2
Lung Cancer 16	Pos	Neg	Neg	Neg	1
Lung Cancer 17	Neg	Neg	Pos	Neg	1
Lung Cancer 12	Neg	Pos	Neg	Neg	1
Lung Cancer 13	Neg	Neg	Pos	Neg	1
Lung Cancer 14	Neg	Neg	Neg	Pos	1
Lung Cancer 17	Neg	Neg	Pos	Neg	
Lung Cancer 18	Neg	Neg	Pos	Neg	
Lung Cancer 20	Neg	Neg	POS	Neg	1
Lung Cancer 22	Neg	Neg	Neg	POS	1
Lung Cancer 20	Pos	Neg	Neg	Neg	1
Apparently Healthy 07	Pos	Pos	Neg	Pos	3
Apparently Healthy 10	Neg	Pos	Neg	Neg	1
Apparently Healthy 16	Neg	Neg	Pos	Neg	1
Apparently Healthy 21	Pos	Neg	Neg	Neg	1
Apparently Healthy 23	Pos	Neg	Neg	Neg	1
Apparently Healthy 38	Neg	Neg	Pos	Neg] 1
Apparently Healthy 45	Neg	Pos	Neg	Neg	1
Apparently Healthy 53	Neg	Neg	Neg	Pos	1
Apparently Healthy 54	Neg	Neg	Neg	Pos	1
Apparently Healthy 56	Neg	Neg	Neg	Pos	1
Apparently Healthy 57	Neg	Neg	Neg	Pos	
Apparently Healthy 77	Pos	Pos	Neg	Neg	2
Apparently Healthy 84	Neg	Neg	POS	Neg	1
Apparently Healthy 88	INEG	Mog	Neg	Neg Nog	
	Neg	Neg	Nog	Doc	1
Δnnarontly Healthy 01	Pos	Pos	Nea	Neg	л Э
	100		ואסע	וזכץ	

Table 2. At the preliminarily selected cutoff for antigen seropositivity of 95% of an apparently healthy population, presumably true positive samples are reactive for several E proteins. In contrast, many of the remaining samples are positive only for one antigen. Larger studies are required to establish criteria to accurately classify true positives, such as requiring positivity for more than one E protein.

Conclusions 6

- preliminarily validated.
- diluted serum or plasma per determination.

7

Acknowledgement

Research reported in this poster was supported by the National Cancer Institute of the National Institutes of Health under Award Number U2CCA271903. The content is solely the responsibility of the authors and does not necessarily represent the official views of the National Institutes of Health.

For Research Use Only. Not for use in diagnostic procedures. $\mathbf{0}$

JRBO-BOOST, TURBO-TAG, N-PLEX, R-PLEX, S-PLEX, T-PLEX, U-PLEX, V-PLEX, MSD (design), MSD (luminous design 1-, 4-, 7-, 9-, & 10-SPOT (designs), 384 WELL 1- & 4-SPOT (designs), N-PLEX (design), R-PLEX (design), U-PLEX (design), V-PLEX (design), V-PLEX (design), It's All About U, Spot the Difference, The Biomarker Company, and The Methodical Mind Experience are service marks owned by or licensed to Meso Scale Diagnostics, LLC. All other trademarks and service marks are the property of their respective owners ©2024 Meso Scale Diagnostics, LLC. All rights reserved

• Sensitive serology assays for oropharyngeal cancer associated HPV-16 E-proteins were developed and

• The assay format is simple, appropriate for high-throughput screening, and uses only 25 µL of 2,500-fold



DOWNLOAD POSTER

