#T2066 Development and Validation of a Human Biomarker 40-Plex V-PLEX[®] Panel – A Tool for Identifying Biomarkers in Health and Disease

Christopher Shelburne, Joseph Manimala, Prachi Gupte, Caitlin Hansen, Amandeep Kaur, Ken Wong, Qian Ning, Allison Cicero, Kristen Weaver, David Stewart, Pankaj Oberoi, and Jacob N. Wohlstadter Meso Scale Discovery, Rockville, Maryland, USA 6 Spike Recovery and Dilution Linearity **1** Abstract **7** Normal Sample Testing **4** Sensitivit

biomarkers that can accurately report biological function or phenotype is necessary for advancing drug concentration corresponding to the signal 2.5 standard deviations above the background (zero calibrator). Limits discovery and diagnostic capabilities. While many biomarker assays are available, their lack of analytical validation and lot-to-lot reproducibility remains a of quantification (LOQ) were determined by spiking a known value of calibrator into diluent to assess the accuracy and precision of the samples across multiple barrier to the success of long-term biomarker programs. These challenges can be addressed with larger multiplex panels that have been validated and lots. The lower LOQ (LLOQ) is the lowest concentration at which the CV of the calculated concentration is <20% and the recovery of the analyte is within 80% to designed to facilitate biomarker discovery in research programs 120% of the known value (75% to 125% for IL-8 in the Chemokine Panel 1 (Human) Kit). The upper LOQ (ULOQ) is the highest concentration at which the CV of Methods: We developed and validated 40 assays for biomarkers relevant to inflammation, immunology, angiogenesis, and vascular injury. These assays the calculated concentration is <20% and the recovery of the analyte is within 80% to 120% of the known value (75% to 125% for IL-8 in the Chemokine Panel 1 detect IFN-γ, IL-1β, IL-2, IL-4, IL-6, IL-8, IL-10, IL-12p70, IL-13, TNFα, GM-CSF, IL-1α, IL-5, IL-7, IL-17A, IL-15, IL-16, IL-12/IL-23p40, TNFβ, IL-8 (high (Human) Kit). The data are from >30 runs collected over 2 months by 4 operators.

abundance), MCP-1, MDC, MCP-4, Eotaxin, IP-10, Eotaxin-3, TARC, MIP-1a, MIP-1B, VEGF, VEGF-C, VEGF-D, Tie-2, sFIt1, PIGF, bFGF, SAA, CRP, VCAM-1 and ICAM-1. All assays were developed with highly characterized and gualified reagents in accordance with fit-for-purpose industry guidelines. Diluents, other kit components, and an optimized assay protocol were developed to help achieve optimum performance. The 40-plex panel rapidly quantified these biomarkers with high precision and accuracy in different human matrices.

Results: We present the assay validation procedure and reproducibility data within and across lots for all 40 assays using a minimum of 3 production lots (generated from independent component lots when possible). The data demonstrate that more than 80% of the assays are sensitive enough to detect low abundance biomarkers in normal biological matrices such as serum, plasma, and urine. Some of the low abundance analytes are quantified at levels below 0.5 pg/ml. In general, these assays display less than 1% non-specific binding with other analytes in the panel or related analytes not in the panel. In addition, we show that subsets of these biomarkers can be multiplexed in any combination without compromising performance

Conclusions: The Human Biomarker 40-Plex V-PLEX Panel provides an accurate and reproducible tool for the identification of a wide array of biomarkers suitable for research or drug development programs. A summary of the development and validation of these panels is presented.

2 Methods

The 40-plex panel was developed as five individual panels designed for optimal performance based on individual assay characteristics. and suitable dilution, diluent, and assay compatibility. The assays presented were developed and validated following fit-for-purpose principles and design control procedures. Accuracy, precision, and specificity were evaluated by multiple operators using independently built kit lots to test calibrators, amples over multiple days and runs. The Pro-inflammatory Panel 1 (human), Cytokine Panel 1 (human), Chemokine Panel 1 (human) Angiogenesis Panel 1 (human), and Vascular Injury Panel 2 (human) kits are based on MSD® ECL technology and follow the same simple three-step



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.

- 1. Add calibrator, control or sample (25 µL/well). Incubate 2 hours at room temperature (RT).
- 2. Wash and add detection antibody solution (25 µL/well). Incubate 1 hour at RT
- 3. Wash and add read buffer (150 µL/well). Analyze with MSD instrument.

3 Calibration Curves

Representative calibration curves are shown below for each of the 40 assays



Pro-inflammatory Panel 1 (Human)			
Analyte	LLOD Median (pg/mL)	LLOQ (pg/mL)	ULOQ (pg/mL)
IFN-γ	0.20	7.47	938
IL-1β	0.04	2.14	375
IL-2	0.09	0.89	938
IL-4	0.02	0.45	158
IL-6	0.06	1.58	488
IL-8	0.04	1.13	375
IL-10	0.03	0.68	233
IL-12p70	0.11	1.22	315
IL-13	0.24	4.21	353
TNFα	0.04	0.69	248

Cytokine Panel 1 (Human)			
Analyte	LLOD Median (pg/mL)	LLOQ (pg/mL)	ULOQ (pg/mL)
GM-CSF	0.14	1.90	750
IL-1α	0.09	2.85	278
IL-5	0.22	6.28	562
IL-7	0.16	1.37	563
IL-12/IL-23p40	0.39	5.68	2250
IL-15	0.17	1.40	525
IL-16	2.83	19.1	1875
IL-17A	0.74	9.32	3653
TNFβ	0.05	1.15	458
VEGF	1.12	7.70	784

Chemokine Panel 1 (Human)			
Analyte	LLOD Median (pg/mL)	LLOQ (pg/mL)	ULOQ (pg/mL)
Eotaxin	3.26	12.3	1120
ΜΙΡ-1β	0.37	2.27	750
Eotaxin-3	1.77	10.2	3750
TARC	0.22	3.32	1120
IP-10	0.37	1.37	500
ΜΙΡ-1α	3.02	13.8	743
IL-8*	95.6	713	43400
MCP-1	0.09	1.09	375
MDC	1.22	88.3	7500
MCP-4	1.69	5.13	469

Because of the high abundance of IL-8 in some sample types Chemokine Panel 1 (Human) uses a low sensitivity II -8 assay

Angiogenesis Panel 1 (Human)			
Analyte	LLOD Median (pg/mL)	LLOQ (pg/mL)	ULOQ (pg/mL)
VEGF-C	11.1	146	17500
VEGF-D	2.53	67.1	18800
Tie-2	21.1	396	63400
Flt-1	0.56	10	6410
PIGF	0.32	10.7	2370
bFGF	0.08	2.6	1780

Vascular Injury Panel 2 (Human)			
Analyte	LLOD Median (pg/mL)	LLOQ (pg/mL)	ULOQ (pg/mL)
SAA	10.9	54.1	137972
CRP	1.33	27.6	49608
VCAM-1	6.00	37.6	31960
ICAM-1	1.03	6.4	42708

5 Specificity and Reproducibility

Each assay in the 40-plex panel was evaluated for cross-reactivity with other analytes in the panel as well as related analytes not in the panel. The cross-reactivity for all assays was <1% (data not shown).



dilution linearity were assessed across multiple lots. Spike recovery testing was conducted in serum, heparin plasma, EDTA plasma and cell culture media. Representative data are presented below showing overall performance of each assay. Almost all assays accurately recovered expected concentrations of spiked analyte in each matrix (dashed lines are at $\pm 20\%$ of expected value).

Spike Recovery











Dilution Linearity

Dilution linearity testing was conducted on all assays. Results were similar with almost all samples recovering within 20% of expected concentrations across all assays. The one exception is PIGF whose performance may benefit from additional dilution.

Normal human matrices (obtained from a commercial vendor) were evaluated for native analyte levels. The representative data for serum is shown with ULOO (dark bars), LLOQ, (black bars) and LLOD (yellow bars) noted for each assay. Nearly all samples were detectable across the 40-plex and many samples were within the LOQ.



8 Whole Blood Testing

Peripheral blood lymphocytes were treated with various stimulants including LPS (10µg/ml), PWM (50µg/ml), anti-CD3/CD28 (5µg/ml each), or ConA (20µg/ml) for 2 hours. The stimulated samples were tested on Pro-inflammatory Panel 1 (human), Cytokine Panel 1 (human) and Chemokine Panel 1 (human) kits. Stimulated studies fo the Angiogenesis Panel 1 (human) and the Vascular Injury Panel 2 (human) were not performed because they are well-represented in normal matrices such as serum Results denoted with (*) are significant relative to controls, and bars denoted with (**) are concentrations above saturation levels. Dose dependent stimulation wa observed for all analytes



9 Conclusion

A 40-analyte panel has been developed and analytically validated for use in the quantification of biologically relevant cytokines and chemokines in multiple human matrices. All assays in these five multiplex panels were individually optimized to achieve the best assay performance and for ease of use, featuring the same simple 3-step protocol requiring just 25 µL of sample to measure all 40 analytes Consistent results were demonstrated for calibration curves and control samples over multiple runs on different days by multiple operators. These assays have sufficient sensitivity, reproducibility, accuracy and performance in different matrices to facilitate biomarker discovery and validation in any research or drug development program.



MESO SCALE DISCOVERY, MESO SCALE DIAGNOSTICS, MSD. DISCOVERY WORKBENCH, MULTI-ARRAY, MULTI-SPOT, QUICKPLEX, SECTOR, SECTOR PR, SECTOR HTS, SULFO-TAG, V-PLE STREPTAVIDIN GOLD, MESO, www.mesoscale.com, SMALL SPOT (design), 96 WELL 1, 4, 7, & 10-SPOT (designs), 384 WELL 1 & 4-SPOT (designs), MSD (design) V-PLEX (design), and SPOT THE DIFFERENCE are trademarks and/or service marks of Meso Scale Diagnostics, LL ©2014 Meso Scale Diagnostics, LLC. All rights reserved.