



CERTIFICATE OF ANALYSIS

Page 1 of 3

COA-05053

Product Description: Proinflammatory Panel 1 (NHP) Kit V-PLEX™ and V-PLEX Plus
Kit Catalog Numbers: K15056D-1, K15056D-2, K15056D-4, K15056G-1, K15056G-2, K15056G-4
Kit Lot Number: K0080259
Expiration Date: 31 AUG 2016

Acceptance Data:

Metric		Specifications	Assay					
			NHP IFN- γ	NHP IL-1 β	NHP IL-2	NHP IL-6	NHP IL-8	NHP IL-10
% Recovery Concentration	Calibrators	80% - 120%	Pass	Pass	Pass	Pass	Pass	Pass
	Controls	80% - 120%	Pass	Pass	Pass	Pass	Pass	Pass
	ULOQ	80% - 120%	Pass	Pass	Pass	Pass	Pass	Pass
	LLOQ	80% - 120%	Pass	Pass	Pass	Pass	Pass	Pass
% Recovery CV	Calibrators	< 20%	Pass	Pass	Pass	Pass	Pass	Pass
	Controls	< 20%	Pass	Pass	Pass	Pass	Pass	Pass
	ULOQ	< 20%	Pass	Pass	Pass	Pass	Pass	Pass
	LLOQ	< 20%	Pass	Pass	Pass	Pass	Pass	Pass
Non-specific binding		< 1%	Pass	Pass	Pass	Pass	Pass	Pass

Concentration of Materials Tested:

Units = pg/mL	Assay					
	NHP IFN- γ *	NHP IL-1 β *	NHP IL-2*	NHP IL-6*	NHP IL-8*	NHP IL-10*
Reconstituted Calibrator	1510	520	1460	749	546	313
Reconstituted Control 1 (A00C0273)	345	129	195	128	129	95.8
Reconstituted Control 2 (A00C0274)	123	52.2	77.6	58.3	37.6	24.7
Reconstituted Control 3 (A00C0293)	37.9	8.62	2.96	8.27	5.19	3.69
ULOQ	1200	415	1160	597	435	250
LLOQ	3.13	1.08	0.890	0.457	0.333	0.191

*The calibrator and controls for this assay contain a human recombinant cytokine calibrator sequence that is highly homologous to the non-human primate cytokine. This assay has been verified to cross-react with homologs of rhesus and cynomolgus monkeys. See Product Insert for experimental data.



CERTIFICATE OF ANALYSIS

Page 2 of 3

COA-05053

Standard Curve Data from Kit QC:

NHP IFN- γ		
Conc., (pg/mL)	Average Signal	%CV
1510	319 790	2.4
378	87 191	2.3
94.4	22 392	2.4
23.6	5683	2.1
5.90	1525	3.6
1.47	543	5.6
0.369	287	10.0
0	215	22.7

NHP IL-1 β		
Conc., (pg/mL)	Average Signal	%CV
520	1 334 727	2.8
130	327 613	3.7
32.5	82 743	2.5
8.13	21 058	4.1
2.03	6311	4.4
0.508	2676	8.8
0.127	1702	4.5
0	1396	9.0

NHP IL-2		
Conc., (pg/mL)	Average Signal	%CV
1460	939 189	2.9
365	261 427	2.8
91.3	69 396	2.8
22.8	16 885	4.0
5.70	4596	3.2
1.43	1385	3.0
0.356	554	4.8
0	284	8.0

NHP IL-6		
Conc., (pg/mL)	Average Signal	%CV
749	1 084 190	3.3
187	246 122	3.3
46.8	55 550	4.1
11.7	12 784	2.4
2.93	3165	3.7
0.731	920	3.4
0.183	343	4.9
0	156	13.0

NHP IL-8		
Conc., (pg/mL)	Average Signal	%CV
546	556 280	2.2
137	141 196	2.5
34.1	35 701	2.3
8.53	8682	2.9
2.13	2346	3.3
0.533	693	5.5
0.133	268	5.9
0	141	13.6

NHP IL-10		
Conc., (pg/mL)	Average Signal	%CV
313	482 307	2.9
78.3	136 010	2.0
19.6	35 672	2.8
4.89	8518	3.6
1.22	2363	2.9
0.306	719	5.4
0.076	323	8.4
0	178	13.7



CERTIFICATE OF ANALYSIS

COA-05053

Proinflammatory Panel 1 (NHP) Kit V-PLEX Components			
Description	Lot Number	Storage Temperature	Expiration Date
Proinflammatory Panel 1 (human) Plate	Z0045508	2-8°C	30 APR 2017
Anti-hu IFN- γ Antibody	D0080363	2-8°C	31 MAR 2017
Anti-hu IL-1 β Antibody	D0080358	2-8°C	30 NOV 2016
Anti-hu IL-2 Antibody	D0080381	2-8°C	31 MAR 2017
Anti-hu IL-6 Antibody	D0080384	2-8°C	30 APR 2017
Anti-hu IL-8 Antibody	D0080380	2-8°C	31 MAR 2017
Anti-hu IL-10 Antibody	D0080298	2-8°C	31 AUG 2016
Proinflammatory Panel 1 (human) Calibrator Blend	A0080045	2-8°C	30 NOV 2019
Diluent 2	M0020085	$\leq -10^{\circ}\text{C}$	31 JAN 2017
Diluent 3	M0030063	$\leq -10^{\circ}\text{C}$	30 APR 2017
MSD Read Buffer T (4X)	Not Kit Specific	Room Temperature	N/A
Additional Components Included in V-PLEX Plus			
Wash Buffer (20X), 100 mL	Not Kit Specific	Room Temperature	N/A
Plate Seals	Not Kit Specific	Room Temperature	N/A
Proinflammatory Panel 1 (human) Control 1	Not Kit Specific; See Separate Control Pack COA		
Proinflammatory Panel 1 (human) Control 2			
Proinflammatory Panel 1 (human) Control 3			

Statement: The above kit lot was tested using MSD procedures and it conforms to MSD’s specifications. The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Caitlin Hansen	Quality Control		19 DEC 2014