



## CERTIFICATE OF ANALYSIS

Page 1 of 2

COA-05092

**Product Description:** Human IP-10 Kit V-PLEX™ and V-PLEX Plus

**Kit Catalog Numbers:** K151NVD-1, K151NVD-2, K151NVD-4, K151NVG-1, K151NVG-2, K151NVG-4

**Kit Lot Number:** K0080264

**Expiration Date:** 31 AUG 2016

### Acceptance Data:

Metric		Specifications	Assay hu IP-10
% Recovery Concentration	Calibrators	80% - 120%	Pass
	Controls	90% - 130%	Pass
	ULOQ	80% - 120%	Pass
	LLOQ	80% - 120%	Pass
% Recovery CV	Calibrators	< 20%	Pass
	Controls	< 20%	Pass
	ULOQ	< 20%	Pass
	LLOQ	< 20%	Pass

### Concentration of Materials Tested:

Units = pg/mL	Assay hu IP-10
Reconstituted Calibrator	2380
Reconstituted Control 1 (A00C0267)	1470
Reconstituted Control 2 (A00C0268)	378
Reconstituted Control 3 (A00C0269)	93.5
ULOQ	500
LLOQ	1.37

### Standard Curve Data from Kit QC:

hu IP-10		
Conc., (pg/mL)	Average Signal	%CV
2380	184 125	8.3
595	102 983	7.1
149	31 777	5.4
37.2	6703	10.3
9.30	2050	3.9
2.32	485	8.7
0.581	165	8.0
0	50	5.6



## CERTIFICATE OF ANALYSIS

Page 2 of 2

COA-05092

Human IP-10 Kit V-PLEX Components			
Description	Lot Number	Storage Temperature	Expiration Date
Human IP-10 Plate	Z0045541	2-8°C	31 MAY 2017
Anti-hu IP-10 Antibody	D0080306	2-8°C	31 AUG 2016
Chemokine Panel 1 (human) Calibrator Blend	A0080037	2-8°C	30 SEP 2016
Diluent 43	M0430011	≤ -10°C	28 FEB 2017
Diluent 3	M0030063	≤ -10°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
Chemokine Panel 1 (human) Control 1	Not Kit Specific; See Separate Control Pack COA		
Chemokine Panel 1 (human) Control 2			
Chemokine 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 JAN 2015