



CERTIFICATE OF ANALYSIS

Page 1 of 2

COA-03790

Product Description: Human MIP-1 β Kit V-PLEX™ and V-PLEX Plus

Kit Catalog Numbers: K151NRD-1, K151NRD-2, K151NRD-4, K151NRG-1, K151NRG-2, K151NRG-4

Kit Lot Number: K0080056

Expiration Date: 30Sep2015

Acceptance Data:

		Assay	
Metric	Specifications	hu MIP-1 β	
% Recovery Concentration	Calibrators	80% - 120%	Pass
	Controls	80% - 120%	Pass
	LLOQ	80% - 120%	Pass
	ULOQ	80% - 120%	Pass
% Recovery CV	Calibrators	<20%	Pass
	Controls	<20%	Pass
	LLOQ	<20%	Pass
	ULOQ	<20%	Pass

Concentration of Materials Tested:

Units = pg/mL	Assay
	hu MIP-1 β
Reconstituted Calibrator	1 100
Reconstituted Control 1 (A00C0190)	678
Reconstituted Control 2 (A00C0191)	175
Reconstituted Control 3 (A00C0192)	36.3
ULOQ	991
LLOQ	2.15

Standard Curve Data from Kit QC:

hu MIP-1 β		
Conc., (pg/ml)	Average Signal	%CV
1 100	356 467	5.2
275	62 404	4.2
68.8	6 689	4.8
17.2	1 035	5.4
4.30	212	6.4
1.07	100	5.1
0.269	70	4.8
0	63	6.1




CERTIFICATE OF ANALYSIS

Page 2 of 2

COA-03790

Human MIP-1 β Kit V-PLEX Components			
Description	Lot Number	Storage Temperature	Expiration Date
Human MIP-1 β Plate	Z0044699	2-8°C	31Oct2015
Anti-hu MIP-1 β Antibody	D0080124	2-8°C	30Dec2015
Chemokine Panel 1 (human) Calibrator Blend	A0080012	2-8°C	30Sep2015
Diluent 43	M0430003	$\leq -10^{\circ}\text{C}$	30Sep2015
Diluent 3	M0030055	$\leq -10^{\circ}\text{C}$	30Sep2015
Read Buffer T (4X)	Not Kit Specific	Room Temperature	N/A
Additional Components Included in V-PLEX Plus			
Wash Buffer (20X), 100mL	Not Kit Specific	Room Temperature	N/A
Plate Seals, pack of 3	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	Paula Eason	Quality Control		29Jul2013