

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Human MIP-3 $\alpha$  Kit  
**Kit Catalog Numbers:** K151XD-Series  
**Kit Lot Number:** K0080620  
**Expiration Date:** 31 Jan 2019

### Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		TH-17 Panel 1 (human) Control 1		TH-17 Panel 1 (human) Control 2		TH-17 Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	hu MIP-3 $\alpha$	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	hu MIP-3 $\alpha$	0.590	80% - 120%	< 20%	325	80% - 120%	< 20%

### Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu MIP-3 $\alpha$	Calibrator	449	pg/mL
		TH-17 Panel 1 (human) Control 1	592	pg/mL
		TH-17 Panel 1 (human) Control 2	235	pg/mL
		TH-17 Panel 1 (human) Control 3	83	pg/mL

Controls tested may not be the same lot as shipped.

### Standard Curve Data:

hu MIP-3 $\alpha$ - Spot 1		
Conc. pg/mL	ECL Signal	%CV
449	177,000	2.4
112	42,700	3.6
28.1	9,660	2.7
7.02	2,320	2.4
1.75	686	2.9
0.438	275	2.8
0.110	178	3.9
0.00	142	3.1



## CERTIFICATE OF ANALYSIS

Page 2 of 2

### Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human MIP-3 $\alpha$ Plate	Z0046183	2-8°C	28 Feb 2019
Anti-hu MIP-3 $\alpha$ Antibody	D0080763	2-8°C	31 Jan 2019
TH17 Panel 1 (human) Calibrator Blend	A0080101	2-8°C	30 Sep 2021
Diluent 43	M0430018	$\leq$ -10°C	31 Jan 2019
Diluent 3	M0030073	$\leq$ -10°C	31 Mar 2019
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
Plate Seals	Not Kit Specific
Wash Buffer (20x), 100mL	Not Kit Specific
TH-17 Panel 1 (human) Control 1	Not Kit Specific
TH-17 Panel 1 (human) Control 2	Not Kit Specific
TH-17 Panel 1 (human) Control 3	Not Kit Specific

### Statement:

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance		03 MAR 2017