

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Human GM-CSF Kit  
**Kit Catalog Numbers:** K156RI-Series, K151RI-Series  
**Kit Lot Number:** K0081504  
**Expiration Date:** 28 Feb 2022

**Acceptance Data/ Percent Recovery Specifications:**

Spot	Name	Calibrator		Cytokine Panel 1 (human) Control 1		Cytokine Panel 1 (human) Control 2		Cytokine Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	GM-CSF	80% - 120%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	GM-CSF	0.842	80% - 120%	< 20%	750	80% - 120%	< 20%

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	GM-CSF	Calibrator	808	pg/mL
		Cytokine Panel 1 (human) Control 1	435	pg/mL
		Cytokine Panel 1 (human) Control 2	69.6	pg/mL
		Cytokine Panel 1 (human) Control 3	11.2	pg/mL

Controls tested may not be the same lot as shipped.

**Standard Curve Data:**

GM-CSF - Spot 1		
Conc. pg/mL	ECL Signal	%CV
808	171,000	1.9
202	45,000	2.5
50.5	11,500	1.9
12.6	2,950	1.6
3.16	782	1.8
0.789	237	2.5
0.197	106	4.3
0.00	66	9.2

## CERTIFICATE OF ANALYSIS


**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human GM-CSF Plate	Z0047350	2-8°C	30 Nov 2022
Anti-hu GM-CSF Antibody	D0081143	2-8°C	28 Feb 2022
Cytokine Panel 1 (human) Calibrator Blend	A0080227	2-8°C	31 Oct 2024
Diluent 43	M0430035	≤ -10°C	31 Mar 2022
Diluent 3	M0030092	≤ -10°C	31 Mar 2022
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
MSD Wash Buffer (20X), 100mL	Not Kit Specific
Plate Seals	Not Kit Specific
Cytokine Panel 1 (human) Control 1	Not Kit Specific
Cytokine Panel 1 (human) Control 2	Not Kit Specific
Cytokine 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>	Karen Hamilla	Quality		20 JUL 2020