

**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:** K0081536  
**Expiration Date:** 30 Jun 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	132,000	4.4
202	33,300	2.7
50.5	8,180	1.1
12.6	2,020	2.7
3.16	548	2.0
0.789	189	3.6
0.197	97	4.4
0.00	67	5.5

**CERTIFICATE OF ANALYSIS**

Page 2 of 2


**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human GM-CSF Plate	Z0047378	2-8°C	28 Feb 2023
Anti-hu GM-CSF Antibody	D0081205	2-8°C	30 Jun 2022
Cytokine Panel 1 (human) Calibrator Blend	A0080227	2-8°C	31 Oct 2024
Diluent 43	M0430037	≤ -10°C	31 Aug 2022
Diluent 3	M0030096	≤ -10°C	30 Jun 2022
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
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		09 OCT 2020