

CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus Human GM-CSF Kit
Kit Catalog Numbers: K151RI-Series, K156RI-Series
Kit Lot Number: K0081289
Expiration Date: 31 Dec 2021

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	1145	pg/mL
		Cytokine Panel 1 (human) Control 1	394	pg/mL
		Cytokine Panel 1 (human) Control 2	61.7	pg/mL
		Cytokine Panel 1 (human) Control 3	8.11	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
1150	146,000	2.1
286	39,500	2.1
71.6	10,300	3.2
17.9	2,620	2.6
4.47	692	4.8
1.12	208	6.2
0.280	96	4.1
0.00	54	8.4

CERTIFICATE OF ANALYSIS

Components:


Description	Lot Number	Storage Temperature	Expiration Date
Human GM-CSF Plate	Z0047141	2-8°C	31 Jan 2022
Anti-hu GM-CSF Antibody	D0081205	2-8°C	30 Jun 2022
Cytokine Panel 1 (human) Calibrator Blend	A0080188	2-8°C	30 Nov 2023
Diluent 43	M0430033	≤ -10°C	31 Jan 2022
Diluent 3	M0030090	≤ -10°C	31 Dec 2021
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
Review/Approval	Karen Hamilla	Quality		13 DEC 2019