

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human IL-10 Kit
Kit Catalog Numbers: K151QU-Series; K156QU-Series
Kit Lot Number: K0081454
Expiration Date: 31 Jul 2022

Acceptance Data/ Percent Recovery Specifications:

		Calibrator		Proinflammatory Panel 1 (human) Control 1		Proinflammatory Panel 1 (human) Control 2		Proinflammatory Panel 1 (human) Control 3	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-10	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	IL-10	0.298	80% - 120%	< 20%	233	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	IL-10	Calibrator	373	pg/mL
		Proinflammatory Panel 1 (human) Control 1	135	pg/mL
		Proinflammatory Panel 1 (human) Control 2	24.8	pg/mL
		Proinflammatory Panel 1 (human) Control 3	3.62	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

IL-10 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
373	246,000	2.7
93.3	64,700	2.7
23.3	16,800	2.8
5.83	4,260	3.7
1.46	1,100	3.6
0.364	329	3.3
0.0911	140	3.0
0.00	77	2.2

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-10 Plate	Z0047294	2-8°C	31 Aug 2022
Anti-hu IL-10 Antibody	D0081437	2-8°C	30 Jun 2023
Proinflammatory Panel 1 (human) Calibrator Blend	A0080216	2-8°C	30 Jun 2024
Diluent 2	M0020112	≤ -10°C	31 Jul 2022
Diluent 3	M0030097	≤ -10°C	31 Aug 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
Proinflammatory Panel 1 (human) Control 1	Not Kit Specific
Proinflammatory Panel 1 (human) Control 2	Not Kit Specific
Proinflammatory 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		26 MAY 2020