

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus® Human IL-6 Kit
Kit Catalog Numbers: K151QX-Series
Kit Lot Number: K0081060
Expiration Date: 31 Jan 2020

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Proinflammatory Panel 1 (human) Control 1		Proinflammatory Panel 1 (human) Control 2		Proinflammatory Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	hu IL-6	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	hu IL-6	0.633	80% - 120%	< 20%	488	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu IL-6	Calibrator	765	pg/mL
		Proinflammatory Panel 1 (human) Control 1	271	pg/mL
		Proinflammatory Panel 1 (human) Control 2	43.5	pg/mL
		Proinflammatory Panel 1 (human) Control 3	8.05	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

hu IL-6 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
765	579,000	3.6
191	136,000	2.5
47.8	29,900	3.7
12.0	7,080	3.4
2.99	1,830	2.2
0.747	504	3.0
0.187	177	2.2
0.00	70	8.2

CERTIFICATE OF ANALYSIS


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-6 Plate	Z0046881	2-8°C	31 Dec 2020
Anti-hu IL-6 Antibody	D0081018	2-8°C	31 Jan 2020
Proinflammatory Panel 1 (human) Calibrator Blend	A0080169	2-8°C	31 May 2023
Diluent 2	M0020099	≤ -70°C	31 Mar 2020
Diluent 3	M0030084	≤ -70°C	31 Jan 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), 100 mL	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	Dawn Stollar	Quality Assurance		24 SEP 2018