

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human IL-6 Kit
Kit Catalog Numbers: K151QX-Series, K156QX-Series
Kit Lot Number: K0081480
Expiration Date: 31 Jul 2022

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	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	IL-6	0.633	80% - 120%	< 20%	488	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	IL-6	Calibrator	697	pg/mL
		Proinflammatory Panel 1 (human) Control 1	289	pg/mL
		Proinflammatory Panel 1 (human) Control 2	49.1	pg/mL
		Proinflammatory Panel 1 (human) Control 3	7.89	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

IL-6 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
697	348,000	1.1
174	80,700	1.1
43.6	18,800	1.4
10.9	4,670	1.9
2.72	1,240	1.8
0.681	356	2.3
0.170	134	5.1
0.00	65	6.3

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-6 Plate	Z0047320	2-8°C	30 Sep 2022
Anti-hu IL-6 Antibody	D0081420	2-8°C	30 Apr 2023
Proinflammatory Panel 1 (human) Calibrator Blend	A0080254	2-8°C	30 Apr 2025
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		28 APR 2020