

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-Plex® and V-Plex Mouse IL-6 Kit
Kit Catalog Numbers: K152QX-Series
Kit Lot Number: K0081242
Expiration Date: 30 Jun 2021

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Proinflammatory Panel 1 (mouse) Control 1		Proinflammatory Panel 1 (mouse) Control 2		Proinflammatory Panel 1 (mouse) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-6	80% - 120%	< 20%	75% - 125%*	< 25%	75% - 125%	< 25%	75% - 125%	< 25%

* See Additional Comments.

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-6	7.61	75% - 125%	< 25%	3140	75% - 125%	< 25%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	IL-6	Calibrator	4730	pg/mL
		Proinflammatory Panel 1 (mouse) Control 1	1050	pg/mL
		Proinflammatory Panel 1 (mouse) Control 2	167	pg/mL
		Proinflammatory Panel 1 (mouse) Control 3	31.5	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
4730	123,000	1.7
1180	26,800	2.5
296	6,470	2.5
73.9	1,580	2.3
18.5	471	2.6
4.62	181	2.8
1.15	113	4.3
0.00	92	8.1

CERTIFICATE OF ANALYSIS

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-6 Plate	Z0047086	2-8°C	30 Sep 2021
Anti-ms IL-6 Antibody	D0081088	2-8°C	31 Oct 2021
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080168	2-8°C	31 May 2023
Diluent 41	M0410027	≤ -10°C	31 Aug 2021
Diluent 45	M0450026	≤ -10°C	30 Jun 2021
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

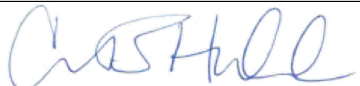
Additional Components Included in V-PLEX Plus	
MSD Wash Buffer (20X), 1000mL	Not Kit Specific
Plate Seals	Not Kit Specific
Proinflammatory Panel 1 (mouse) Control 1	Not Kit Specific
Proinflammatory Panel 1 (mouse) Control 2	Not Kit Specific
Proinflammatory Panel 1 (mouse) Control 3	Not Kit Specific

Additional Comments:

	Assay	Spot	Specification	Comment
1	IL-6	1	Control Conc.	Control-1 recovery 129.4%

Statement:

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Carissa Hubbard	Quality		02 JUL 2019