

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Mouse IL-10 Kit
Kit Catalog Numbers: K152QU-series
Kit Lot Number: K0080790
Expiration Date: 31 May 2019

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	ms IL-10	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	ms IL-10	19.8	75% - 125%	< 25%	2030	75% - 125%	< 25%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	ms IL-10	Calibrator	3070	pg/mL
		Proinflammatory Panel 1 (mouse) Control 1	1040	pg/mL
		Proinflammatory Panel 1 (mouse) Control 2	312	pg/mL
		Proinflammatory Panel 1 (mouse) Control 3	97.7	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

ms IL-10 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
3070	489,000	2.8
768	114,000	2.5
192	28,100	2.3
48.0	6,810	4.0
12.0	1,810	3.0
3.00	541	2.1
0.750	228	4.8
0.00	120	6.9

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-10 Plate	Z0046510	2-8°C	31 Oct 2019
Anti-ms IL-10 Antibody	D0080858	2-8°C	31 May 2019
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080110	2-8°C	31 Dec 2021
Diluent 41	M0410017	≤ -10°C	31 Oct 2019
Diluent 45	M0450018	≤ -10°C	30 Nov 2019
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 (mouse) Control 1	Not Kit Specific
Proinflammatory Panel 1 (mouse) Control 2	Not Kit Specific
Proinflammatory Panel 1 (mouse) 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	Vanessa Castleberry	Quality Assurance	<i>Vanessa Castleberry</i>	28AUG2017