

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Mouse IL-6 Kit
Kit Catalog Numbers: K152QX-series
Kit Lot Number: K0080817
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-6	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-6	7.61	75% - 125%	< 25%	3140	75% - 125%	< 25%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	ms IL-6	Calibrator	5200	pg/mL
		Proinflammatory Panel 1 (mouse) Control 1	1590	pg/mL
		Proinflammatory Panel 1 (mouse) Control 2	247	pg/mL
		Proinflammatory Panel 1 (mouse) Control 3	40.1	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

ms IL-6 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
5200	225,000	2.7
1300	53,100	2.0
325	13,000	2.8
81.3	3,310	3.2
20.3	948	4.2
5.08	305	2.8
1.27	143	3.5
0.00	99	7.3

CERTIFICATE OF ANALYSIS

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-6 Plate	Z0046518	2-8°C	31 Oct 2019
Anti-ms IL-6 Antibody	D0080856	2-8°C	31 May 2019
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080060	2-8°C	31 Jul 2020
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