

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human IL-2 Kit
Kit Catalog Numbers: K151QQ-series
Kit Lot Number: K0080865
Expiration Date: 31 Oct 2019

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-2	80% - 120%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	hu IL-2	0.890	80% - 120%	< 20%	938	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu IL-2	Calibrator	1520	pg/mL
		Proinflammatory Panel 1 (human) Control 1	431	pg/mL
		Proinflammatory Panel 1 (human) Control 2	42.3	pg/mL
		Proinflammatory Panel 1 (human) Control 3	3.83	pg/mL

Controls tested may not be the same lot as shipped.

CERTIFICATE OF ANALYSIS

Standard Curve Data:

hu IL-2 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1520	499,000	1.6
380	138,000	1.8
95.0	36,900	1.5
23.8	9,280	3.4
5.94	2,370	3.2
1.48	635	2.0
0.371	206	3.9
0.00	65	4.7

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-2 Plate	Z0046587	2-8°C	31 Dec 2019
Anti-hu IL-2 Antibody	D0080982	2-8°C	31 Oct 2019
Proinflammatory Panel 1 (human) Calibrator Blend	A0080139	2-8°C	31 Jul 2022
Diluent 2	M0020096	≤ -10°C	30 Nov 2019
Diluent 3	M0030077	≤ -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 (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	Vanessa Castleberry	Quality Assurance	<i>Vanessa Castleberry</i>	29AUG2017