

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Mouse IL-27p28/IL-30 Kit
Kit Catalog Numbers: K152WX-Series
Kit Lot Number: K0080725
Expiration Date: 31 Aug 2019

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Cytokine Panel (mouse) Control 1		Cytokine Panel (mouse) Control 2		Cytokine Panel (mouse) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	ms IL-27p28/IL-30	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	ms IL-27p28/IL-30	5.91	75% - 125%	< 20%	6500	75% - 125%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	ms IL-27p28/IL-30	Calibrator	10500	pg/mL
		Cytokine Panel (mouse) Control 1	8930	pg/mL
		Cytokine Panel (mouse) Control 2	1290	pg/mL
		Cytokine Panel (mouse) Control 3	203	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

ms IL-27p28/IL-30 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
10500	415,000	2.5
2630	126,000	3.5
656	31,600	3.3
164	7,510	2.0
41.0	1,940	2.3
10.3	580	1.5
2.56	244	3.4
0.00	119	4.2

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-27p28/IL-30 Plate	Z0046392	2-8°C	31 Aug 2019
Anti-ms IL-27p28/IL-30 Antibody	D0080937	2-8°C	31 Aug 2019
Cytokine Panel 1 (mouse) Calibrator Blend	A0080125	2-8°C	31 Mar 2022
Diluent 41	M0410016	≤ -10°C	31 Aug 2019
Diluent 45	M0450016	≤ -10°C	30 Sep 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
Cytokine Panel (mouse) Control 1	Not Kit Specific
Cytokine Panel (mouse) Control 2	Not Kit Specific
Cytokine Panel (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	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	23 JUN 2017