



RFE MEMO

Page 1 of 1

Product Description(s): V-PLEX® and V-PLEX Plus Cytokine Panel 2 (human) Kit

Catalog Number(s) / Size(s): K15084-Series

Lot Number: K0080804

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0080804 has been extended to 30 APR 2020.

The expiration dates for the following kit components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Cytokine Panel 2 (human) Plate	Z0046488	31 OCT 2019	30 APR 2020
Anti-hu IL-17 A/F Antibody	D0080766	30 APR 2019	30 APR 2020
Anti-hu IL-17B Antibody	D0080767	30 APR 2019	30 APR 2020
Anti-hu IL-17C Antibody	D0080768	30 APR 2019	30 APR 2020
Anti-hu IL-17D Antibody	D0080946	31 AUG 2019	30 APR 2020
Anti-hu IL-IRA Antibody	D0080771	30 APR 2019	30 APR 2020
Anti-hu IL-3 Antibody	D0080772	30 APR 2019	30 APR 2020
Anti-hu IL-9 Antibody	D0080773	30 APR 2019	30 APR 2020
Anti-hu TSLP Antibody	D0080775	30 APR 2019	30 APR 2020
Diluent 43	M0430021	30 SEP 2019	30 APR 2020
Diluent 3	M0030077	30 NOV 2019	30 APR 2020

Function	Name	Signature	Date
Quality Assurance	Dawn Stollar	<i>Dawn Stollar</i>	29 JAN 2019

This memo supersedes all other statements regarding date of expiration for this product.

CERTIFICATE OF ANALYSIS

Page 1 of 4

Product Description: V-PLEX® and V-PLEX Plus Cytokine Panel 2 (human) Kit

Kit Catalog Numbers: K15084-Series

Kit Lot Number: K0080804

Expiration Date: 30 Apr 2019

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Cytokine Panel 2 (human) Control 1		Cytokine Panel 2 (human) Control 2		Cytokine Panel 2 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	hu IL-17 A/F	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
2	hu IL-17B	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
3	hu IL-17C	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
4	hu IL-17D	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
6	hu IL-1RA	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
7	hu IL-3	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
8	hu IL-9	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
10	hu TSLP	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	hu IL-17 A/F	7.57	80% - 120%	< 20%	3900	80% - 120%	< 20%
2	hu IL-17B	1.12	75% - 125%	< 20%	1040	75% - 125%	< 20%
3	hu IL-17C	3.84	75% - 125%	< 20%	1620	75% - 125%	< 20%
4	hu IL-17D	11.2	75% - 125%	< 20%	5200	75% - 125%	< 20%
6	hu IL-1RA	9.19	75% - 125%	< 25%	650	80% - 120%	< 20%
7	hu IL-3	12.6	75% - 125%	< 20%	1950	80% - 120%	< 20%
8	hu IL-9	2.23	80% - 120%	< 25%	975	80% - 120%	< 20%
10	hu TSLP	0.460	75% - 125%	< 20%	325	75% - 125%	< 20%

Acceptance Data/ Non Specific Binding Specifications:

Spot	Name	NSB Spec
1	hu IL-17 A/F	< 1.0%
2	hu IL-17B	< 1.0%
3	hu IL-17C	< 1.0%
4	hu IL-17D	< 1.0%
6	hu IL-1RA	< 1.0%
7	hu IL-3	< 1.0%
8	hu IL-9	< 1.0%
10	hu TSLP	< 1.0%

CERTIFICATE OF ANALYSIS

Page 2 of 4

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu IL-17 A/F	Calibrator	5607	pg/mL
		Cytokine Panel 2 (human) Control 1	5913	pg/mL
		Cytokine Panel 2 (human) Control 2	796	pg/mL
		Cytokine Panel 2 (human) Control 3	74	pg/mL
2	hu IL-17B	Calibrator	1520	pg/mL
		Cytokine Panel 2 (human) Control 1	487	pg/mL
		Cytokine Panel 2 (human) Control 2	71	pg/mL
		Cytokine Panel 2 (human) Control 3	133	pg/mL
3	hu IL-17C	Calibrator	2627	pg/mL
		Cytokine Panel 2 (human) Control 1	2360	pg/mL
		Cytokine Panel 2 (human) Control 2	295	pg/mL
		Cytokine Panel 2 (human) Control 3	77	pg/mL
4	hu IL-17D	Calibrator	7467	pg/mL
		Cytokine Panel 2 (human) Control 1	5479	pg/mL
		Cytokine Panel 2 (human) Control 2	559	pg/mL
		Cytokine Panel 2 (human) Control 3	146	pg/mL
6	hu IL-1RA	Calibrator	732	pg/mL
		Cytokine Panel 2 (human) Control 1	670	pg/mL
		Cytokine Panel 2 (human) Control 2	143	pg/mL
		Cytokine Panel 2 (human) Control 3	50	pg/mL
7	hu IL-3	Calibrator	3060	pg/mL
		Cytokine Panel 2 (human) Control 1	3173	pg/mL
		Cytokine Panel 2 (human) Control 2	703	pg/mL
		Cytokine Panel 2 (human) Control 3	240	pg/mL
8	hu IL-9	Calibrator	1490	pg/mL
		Cytokine Panel 2 (human) Control 1	1272	pg/mL
		Cytokine Panel 2 (human) Control 2	345	pg/mL
		Cytokine Panel 2 (human) Control 3	66	pg/mL
10	hu TSLP	Calibrator	563	pg/mL
		Cytokine Panel 2 (human) Control 1	581	pg/mL
		Cytokine Panel 2 (human) Control 2	87	pg/mL
		Cytokine Panel 2 (human) Control 3	12	pg/mL

Controls tested may not be the same lot as shipped.

CERTIFICATE OF ANALYSIS

Page 3 of 4

Standard Curve Data:

hu IL-17 A/F - Spot 1			hu IL-17B - Spot 2			hu IL-17C - Spot 3		
Conc. pg/mL	ECL Signal	%CV	Conc. pg/mL	ECL Signal	%CV	Conc. pg/mL	ECL Signal	%CV
5610	835,000	3.0	1520	830,000	2.5	2630	326,000	8.2
1400	214,000	2.3	380	209,000	2.9	657	85,400	7.5
350	50,900	2.9	95.0	51,000	3.3	164	22,300	7.1
87.6	12,300	1.9	23.8	12,200	2.7	41.0	5,420	6.3
21.9	3,390	1.8	5.94	3,180	2.4	10.3	1,570	6.6
5.48	1,110	3.4	1.48	934	3.1	2.57	567	4.3
1.37	499	7.2	0.371	337	7.5	0.641	312	9.3
0.00	293	14.0	0.00	134	18.0	0.00	215	14.0

hu IL-17D - Spot 4			hu IL-1RA - Spot 6			hu IL-3 - Spot 7		
Conc. pg/mL	ECL Signal	%CV	Conc. pg/mL	ECL Signal	%CV	Conc. pg/mL	ECL Signal	%CV
7470	289,000	7.7	732	576,000	4.1	3060	956,000	3.8
1870	82,000	5.8	183	109,000	3.8	765	180,000	2.5
467	21,000	5.9	45.8	20,700	4.0	191	24,300	3.8
117	5,320	4.1	11.4	4,660	5.4	47.8	3,570	3.8
29.2	1,520	4.0	2.86	1,560	7.8	12.0	908	4.1
7.29	587	5.0	0.715	935	21.0	2.99	353	10.0
1.82	322	8.7	0.179	786	11.0	0.747	233	17.0
0.00	259	16.0	0.00	556	25.0	0.00	189	17.0

hu IL-9 - Spot 8			hu TSLP - Spot 10		
Conc. pg/mL	ECL Signal	%CV	Conc. pg/mL	ECL Signal	%CV
1490	1,330,000	5.2	563	666,000	2.7
373	340,000	4.5	141	195,000	2.3
93.1	86,700	4.9	35.2	51,700	3.2
23.3	21,800	4.1	8.80	12,500	4.0
5.82	5,640	5.3	2.20	3,320	2.7
1.46	1,660	5.8	0.550	978	5.0
0.364	615	7.9	0.137	378	8.3
0.00	273	30.0	0.00	178	17.0

CERTIFICATE OF ANALYSIS

Page 4 of 4

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Cytokine Panel 2 (human) Plate	Z0046488	2-8°C	31 Oct 2019
Anti-hu IL-17 A/F Antibody	D0080766	2-8°C	30 Apr 2019
Anti-hu IL-17B Antibody	D0080767	2-8°C	30 Apr 2019
Anti-hu IL-17C Antibody	D0080768	2-8°C	30 Apr 2019
Anti-hu IL-17D Antibody	D0080946	2-8°C	31 Aug 2019
Anti-hu IL-1RA Antibody	D0080771	2-8°C	30 Apr 2019
Anti-hu IL-3 Antibody	D0080772	2-8°C	30 Apr 2019
Anti-hu IL-9 antibody	D0080773	2-8°C	30 Apr 2019
Anti-hu TSLP Antibody	D0080775	2-8°C	30 Apr 2019
Cytokine Panel 2 (human) Calibrator Blend	A0080126	2-8°C	31 Mar 2022
Diluent 43	M0430021	≤ -10°C	30 Sep 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
Cytokine Panel 2 (human) Control 1	Not Kit Specific
Cytokine Panel 2 (human) Control 2	Not Kit Specific
Cytokine Panel 2 (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	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	27 JUL 2017