



RFE MEMO

Page 1 of 1

Product Description(s): V-PLEX® and V-PLEX Plus TH17 Panel 1(mouse) Kit

Catalog Number(s) / Size(s): K15246-Series, K152WU-Series

Lot Number: K0081130

Based on additional testing, the original date of expiration for this kit has been extended by 12 months. The expiration date of K0081130 has been extended to 31 OCT 2020.

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

Description	Lot Number	Original Expiration Date	New Expiration Date
Anti-ms MIP-3 α Antibody	D0081048	31 JUL 2020	31 OCT 2020
Anti-ms IL-23 Antibody	D0080951	30 NOV 2019	30 NOV 2020
Anti-ms IL-17C Antibody	D0080952	30 NOV 2019	30 NOV 2020
Anti-ms IL-31 Antibody	D0080953	30 NOV 2019	30 NOV 2020
Anti-ms IL-21 Antibody	D0080954	31 OCT 2019	31 OCT 2020
Anti-ms IL-16 Antibody	D0081050	30 APR 2020	31 OCT 2020
Anti-ms IL-17E/IL-25 Antibody	D0081051	30 JUN 2020	31 OCT 2020

Function	Name	Signature	Date
Quality	Xiomara Talavera		20 SEP 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 TH17 Panel 1(mouse) Kit

Kit Catalog Numbers: K15246-Series, K152WU-Series

Kit Lot Number: K0081130

Expiration Date: 31 Oct 2019

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		TH17 Panel 1(mouse) Control 1		TH17 Panel 1(mouse) Control 2		TH17 Panel 1(mouse) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	ms MIP-3α	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
2	ms IL-22	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
3	ms IL-23	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
4	ms IL-17C	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
5	ms IL-31	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
6	ms IL-21	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
8	ms IL-16	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
9	ms IL-17A	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
10	ms IL-17E/IL-25	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	ms MIP-3α	3.42	75% - 125%	< 20%	530	75% - 125%	< 20%
2	ms IL-22	1.58	75% - 125%	< 20%	380	75% - 125%	< 20%
3	ms IL-23	4.19	75% - 125%	< 25%	7600	75% - 125%	< 20%
4	ms IL-17C	0.971	75% - 125%	< 20%	1200	75% - 125%	< 20%
5	ms IL-31	22.6	75% - 125%	< 20%	21000	75% - 125%	< 20%
6	ms IL-21	12.1	75% - 125%	< 25%	9300	75% - 125%	< 20%
8	ms IL-16	15.0	75% - 125%	< 25%	1800	75% - 125%	< 20%
9	ms IL-17A	0.255	75% - 125%	< 20%	360	75% - 125%	< 20%
10	ms IL-17E/IL-25	4.54	75% - 125%	< 20%	3300	75% - 125%	< 20%

Acceptance Data/ Non Specific Binding Specifications:

Spot	Name	NSB Spec
1	ms MIP-3α	< 1.0%
2	ms IL-22	< 1.0%
3	ms IL-23	< 1.0%
4	ms IL-17C	< 1.0%
5	ms IL-31	< 1.0%
6	ms IL-21	< 1.0%
8	ms IL-16	< 1.0%
9	ms IL-17A	< 1.0%
10	ms IL-17E/IL-25	< 1.0%

CERTIFICATE OF ANALYSIS

Page 2 of 4

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	ms MIP-3 α	Calibrator	1010	pg/mL
		TH17 Panel 1(mouse) Control 1	797	pg/mL
		TH17 Panel 1(mouse) Control 2	200	pg/mL
		TH17 Panel 1(mouse) Control 3	54.6	pg/mL
2	ms IL-22	Calibrator	573	pg/mL
		TH17 Panel 1(mouse) Control 1	423	pg/mL
		TH17 Panel 1(mouse) Control 2	133	pg/mL
		TH17 Panel 1(mouse) Control 3	23.2	pg/mL
3	ms IL-23	Calibrator	8650	pg/mL
		TH17 Panel 1(mouse) Control 1	6450	pg/mL
		TH17 Panel 1(mouse) Control 2	1330	pg/mL
		TH17 Panel 1(mouse) Control 3	199	pg/mL
4	ms IL-17C	Calibrator	1390	pg/mL
		TH17 Panel 1(mouse) Control 1	1300	pg/mL
		TH17 Panel 1(mouse) Control 2	203	pg/mL
		TH17 Panel 1(mouse) Control 3	35.8	pg/mL
5	ms IL-31	Calibrator	32700	pg/mL
		TH17 Panel 1(mouse) Control 1	24500	pg/mL
		TH17 Panel 1(mouse) Control 2	4990	pg/mL
		TH17 Panel 1(mouse) Control 3	1140	pg/mL
6	ms IL-21	Calibrator	15400	pg/mL
		TH17 Panel 1(mouse) Control 1	11200	pg/mL
		TH17 Panel 1(mouse) Control 2	2100	pg/mL
		TH17 Panel 1(mouse) Control 3	355	pg/mL
8	ms IL-16	Calibrator	2070	pg/mL
		TH17 Panel 1(mouse) Control 1	1810	pg/mL
		TH17 Panel 1(mouse) Control 2	1010	pg/mL
		TH17 Panel 1(mouse) Control 3	548	pg/mL
9	ms IL-17A	Calibrator	638	pg/mL
		TH17 Panel 1(mouse) Control 1	508	pg/mL
		TH17 Panel 1(mouse) Control 2	85.2	pg/mL
		TH17 Panel 1(mouse) Control 3	14.8	pg/mL
10	ms IL-17E/IL-25	Calibrator	5150	pg/mL
		TH17 Panel 1(mouse) Control 1	4230	pg/mL
		TH17 Panel 1(mouse) Control 2	1060	pg/mL
		TH17 Panel 1(mouse) Control 3	283	pg/mL

Controls tested may not be the same lot as shipped.

CERTIFICATE OF ANALYSIS

Standard Curve Data:

ms MIP-3a - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1010	573,000	4.6
253	103,000	5.0
63.1	18,800	5.3
15.8	4,900	4.3
3.95	1,400	4.6
0.986	559	5.8
0.247	328	11.0
0.00	256	8.1

ms IL-22 - Spot 2		
Conc. pg/mL	ECL Signal	%CV
573	1,330,000	2.1
143	282,000	2.0
35.8	56,800	3.9
8.95	13,000	2.3
2.24	3,600	5.5
0.560	1,040	2.4
0.140	456	4.7
0.00	251	3.7

ms IL-23 - Spot 3		
Conc. pg/mL	ECL Signal	%CV
8650	792,000	2.3
2160	273,000	1.7
541	73,200	3.8
135	18,900	2.9
33.8	5,080	2.5
8.45	1,430	3.1
2.11	514	5.7
0.00	216	7.8

ms IL-17C - Spot 4		
Conc. pg/mL	ECL Signal	%CV
1390	524,000	3.6
348	123,000	2.9
86.9	29,500	3.1
21.7	7,530	3.0
5.43	2,290	3.4
1.36	915	2.8
0.339	557	4.0
0.00	434	7.8

ms IL-31 - Spot 5		
Conc. pg/mL	ECL Signal	%CV
32700	863,000	3.5
8180	315,000	2.0
2040	64,500	2.7
511	13,000	2.9
128	3,630	2.1
31.9	1,170	2.4
7.98	592	4.0
0.00	396	4.6

ms IL-21 - Spot 6		
Conc. pg/mL	ECL Signal	%CV
15400	391,000	2.4
3850	128,000	4.5
963	32,800	4.9
241	8,610	5.1
60.2	2,480	3.7
15.0	901	3.3
3.76	495	4.6
0.00	356	4.9

ms IL-16 - Spot 8		
Conc. pg/mL	ECL Signal	%CV
2070	467,000	5.2
518	83,300	3.0
129	14,400	6.0
32.3	3,210	7.9
8.09	1,530	9.1
2.02	1,060	14.0
0.505	952	15.0
0.00	872	14.0

ms IL-17A - Spot 9		
Conc. pg/mL	ECL Signal	%CV
638	841,000	1.9
160	236,000	1.4
39.9	59,600	2.9
9.97	14,900	2.7
2.49	3,990	2.3
0.623	1,150	2.6
0.156	456	5.3
0.00	225	7.0

ms IL-17E/IL-25 - Spot 10		
Conc. pg/mL	ECL Signal	%CV
5150	1,290,000	3.0
1290	275,000	2.2
322	57,000	6.1
80.5	14,300	7.2
20.1	4,000	3.1
5.03	1,300	5.6
1.26	548	5.4
0.00	326	5.4

CERTIFICATE OF ANALYSIS

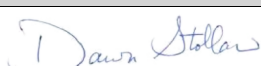
Components:

Description	Lot Number	Storage Temperature	Expiration Date
TH17 Panel 1(mouse) Plate	Z0046952	2-8°C	30 Apr 2021
Anti-ms MIP-3α Antibody	D0081048	2-8°C	31 Jul 2020
Anti-ms IL-22 Antibody	D0081100	2-8°C	31 Aug 2022
Anti-ms IL-23 Antibody	D0080951	2-8°C	30 Nov 2019
Anti-ms IL-17C Antibody	D0080952	2-8°C	30 Nov 2019
Anti-ms IL-31 Antibody	D0080953	2-8°C	30 Nov 2019
Anti-ms IL-21 Antibody	D0080954	2-8°C	31 Oct 2019
Anti-ms IL-16 Antibody	D0081050	2-8°C	30 Apr 2020
Anti-ms IL-17A Antibody	D0081190	2-8°C	31 Jul 2022
Anti-ms IL-17E/IL-25 Antibody	D0081051	2-8°C	30 Jun 2020
TH17 Panel 1 (mouse) Calibrator Blend	A0080157	2-8°C	31 Oct 2022
Diluent 41	M0410025	≤ -10°C	31 Jan 2021
Diluent 45	M0450023	≤ -10°C	31 Jan 2021
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
Wash Buffer (20X), 100mL	Not Kit Specific
Plate Seals	Not Kit Specific
TH17 Panel 1(mouse) Control 1	Not Kit Specific
TH17 Panel 1(mouse) Control 2	Not Kit Specific
TH17 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	Dawn Stollar	Quality Assurance		08 APR 2019