

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human TSLP Kit

Kit Catalog Numbers: K151XE-Series

Kit Lot Number: K0081247

Expiration Date: 31 Dec 2020

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	TSLP	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	TSLP	0.460	75% - 125%	< 20%	325	75% - 125%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	TSLP	Calibrator	515	pg/mL
		Cytokine Panel 2 (human) Control 1	619	pg/mL
		Cytokine Panel 2 (human) Control 2	71.5	pg/mL
		Cytokine Panel 2 (human) Control 3	12.1	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

TSLP - Spot 1		
Conc. pg/mL	ECL Signal	%CV
515	180,000	3.1
129	31,900	6.4
32.2	7,200	9.2
8.05	1,780	7.6
2.01	505	6.0
0.503	181	3.4
0.126	110	3.1
0.00	85	7.5

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human TSLP Plate	Z0047093	2-8°C	31 Oct 2021
Anti-hu TSLP Antibody	D0081301	2-8°C	31 Dec 2022
Cytokine Panel 2 (human) Calibrator Blend	A0080116	2-8°C	31 Jan 2022
Diluent 43	M0430027	≤ -10°C	31 Dec 2020
Diluent 3	M0030084	≤ -10°C	31 Jan 2021
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), 100 mL	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	Karen Hamilla	Quality		28 OCT 2019