

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human Eotaxin Kit
Kit Catalog Numbers: K151NS-Series
Kit Lot Number: K0080910
Expiration Date: 31 Jul 2019

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Chemokine Panel 1 (human) Control 1		Chemokine Panel 1 (human) Control 2		Chemokine Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	hu Eotaxin	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	hu Eotaxin	12.3	80% - 120%	< 20%	1120	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu Eotaxin	Calibrator	1690	pg/mL
		Chemokine Panel 1 (human) Control 1	1370	pg/mL
		Chemokine Panel 1 (human) Control 2	518	pg/mL
		Chemokine Panel 1 (human) Control 3	160	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

hu Eotaxin - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1690	339,000	1.0
423	63,600	2.2
106	9,080	1.6
26.4	1,340	2.0
6.60	501	7.0
1.65	168	3.7
0.413	90	4.4
0.00	66	3.9



CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human Eotaxin Plate	Z0046684	2-8°C	30 Apr 2020
Anti-hu Eotaxin Antibody	D0080897	2-8°C	31 Jul 2019
Chemokine Panel 1 (human) Calibrator Blend	A0080079	2-8°C	31 May 2021
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
Chemokine Panel 1 (human) Control 1	Not Kit Specific
Chemokine Panel 1 (human) Control 2	Not Kit Specific
Chemokine 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	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	27 NOV 2017