

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human Eotaxin-3 Kit
Kit Catalog Numbers: K151NU-series
Kit Lot Number: K0080714
Expiration Date: 30 Jun 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-3	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-3	10.2	80% - 120%	< 20%	3750	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu Eotaxin-3	Calibrator	7520	pg/mL
		Chemokine Panel 1 (human) Control 1	4670	pg/mL
		Chemokine Panel 1 (human) Control 2	794	pg/mL
		Chemokine Panel 1 (human) Control 3	127	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

hu Eotaxin-3 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
7520	61,200	3.2
1880	20,600	3.6
470	6,070	4.2
118	1,640	1.4
29.4	456	2.6
7.34	158	4.0
1.84	87	4.0
0.00	61	5.0

CERTIFICATE OF ANALYSIS

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human Eotaxin-3 Plate	Z0046376	2-8°C	31 Jul 2019
Anti-hu Eotaxin-3 Antibody	D0080862	2-8°C	30 Jun 2019
Chemokine Panel 1 (human) Calibrator Blend	A0080106	2-8°C	30 Nov 2021
Diluent 43	M0430020	≤ -10°C	30 Jun 2019
Diluent 3	M0030074	≤ -10°C	30 Jun 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>	28 FEB 2017