

**CERTIFICATE OF ANALYSIS**

**Product Description:** V-PLEX® and V-PLEX Plus Human Eotaxin-3 Kit  
**Kit Catalog Numbers:** K151NU-series  
**Kit Lot Number:** K0080728  
**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	7120	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.

**CERTIFICATE OF ANALYSIS**

Page 2 of 2

**Standard Curve Data:**

hu Eotaxin-3 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
7120	69,000	3.1
1780	21,500	3.0
445	5,850	3.1
111	1,530	1.8
27.8	429	1.3
6.95	149	3.3
1.74	78	1.5
0.00	57	5.4


**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human Eotaxin-3 Plate	Z0046404	2-8°C	31 Jul 2019
Anti-hu Eotaxin-3 Antibody	D0080862	2-8°C	30 Jun 2019
Chemokine Panel 1 (human) Calibrator Blend	A0080122	2-8°C	28 Feb 2022
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
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance		27 MAR 2017