

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human Eotaxin-3 Kit

Kit Catalog Numbers: K151NU-Series, K156NU-Series

Kit Lot Number: K0081533

Expiration Date: 31 Aug 2022

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	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	Eotaxin-3	10.2	80% - 120%	< 20%	3750	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	Eotaxin-3	Calibrator	4990	pg/mL
		Chemokine Panel 1 (human) Control 1	4250	pg/mL
		Chemokine Panel 1 (human) Control 2	593	pg/mL
		Chemokine Panel 1 (human) Control 3	94.6	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

Eotaxin-3 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
4990	90,500	4.5
1250	23,200	3.4
312	6,260	2.0
78.0	1,610	2.2
19.5	448	2.7
4.87	155	3.3
1.22	84	5.2
0.00	61	5.9

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human Eotaxin-3 Plate	Z0047372	2-8°C	31 Jan 2023
Anti-hu Eotaxin-3 Antibody	D0081260	2-8°C	31 Oct 2022
Chemokine Panel 1 (human) Calibrator Blend	A0080259	2-8°C	31 May 2025
Diluent 43	M0430039	≤ -10°C	30 Sep 2022
Diluent 3	M0030097	≤ -10°C	31 Aug 2022
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	Karen Hamilla	Quality		18 SEP 2020