

**CERTIFICATE OF ANALYSIS**

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Human IL-1 $\beta$  Kit

**Kit Catalog Numbers:** K151QP-Series, K156QP-Series

**Kit Lot Number:** K0081353

**Expiration Date:** 31 Dec 2020

**Acceptance Data/ Percent Recovery Specifications:**

Spot	Name	Calibrator		Proinflammatory Panel 1 (human) Control 1		Proinflammatory Panel 1 (human) Control 2		Proinflammatory Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-1 $\beta$	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-1 $\beta$	2.14	80% - 120%	< 20%	375	80% - 120%	< 20%

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	IL-1 $\beta$	Calibrator	587	pg/mL
		Proinflammatory Panel 1 (human) Control 1	213	pg/mL
		Proinflammatory Panel 1 (human) Control 2	52	pg/mL
		Proinflammatory Panel 1 (human) Control 3	10.8	pg/mL

Controls tested may not be the same lot as shipped.

**Standard Curve Data:**

IL-1 $\beta$ - Spot 1		
Conc. pg/mL	ECL Signal	%CV
587	607,000	2.4
147	150,000	2.2
36.7	37,000	1.5
9.17	9,050	1.1
2.29	2,380	1.6
0.573	645	2.4
0.143	235	4.0
0.00	98	8.1

## CERTIFICATE OF ANALYSIS

Page 2 of 2


**Components:**

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-1 $\beta$ Plate	Z0046863	2-8°C	31 Dec 2020
Anti-hu IL-1 $\beta$ Antibody	D0081082	2-8°C	30 Sep 2021
Proinflammatory Panel 1 (human) Calibrator Blend	A0080153	2-8°C	31 Oct 2022
Diluent 2	M0020108	$\leq$ -10°C	28 Feb 2022
Diluent 3	M0030090	$\leq$ -10°C	31 Dec 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), 100mL	Not Kit Specific
Proinflammatory Panel 1 (human) Control 1	Not Kit Specific
Proinflammatory Panel 1 (human) Control 2	Not Kit Specific
Proinflammatory 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>	Karen Hamilla	Quality		12 FEB 2020