

# CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus Mouse IL-1 $\beta$  Kit  
**Kit Catalog Numbers:** K152QP-Series  
**Kit Lot Number:** K0081239  
**Expiration Date:** 30 Jun 2021

**Acceptance Data/ Percent Recovery Specifications:**

Spot	Name	Calibrator		Proinflammatory Panel 1 (mouse) Control 1		Proinflammatory Panel 1 (mouse) Control 2		Proinflammatory Panel 1 (mouse) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-1 $\beta$	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-1 $\beta$	0.720	75% - 125%	< 25%	1030	75% - 125%	< 25%

**Stock Concentration of Material Tested :**

Spot	Name	Item	Conc	Units
1	IL-1 $\beta$	Calibrator	1430	pg/mL
		Proinflammatory Panel 1 (mouse) Control 1	741	pg/mL
		Proinflammatory Panel 1 (mouse) Control 2	53.8	pg/mL
		Proinflammatory Panel 1 (mouse) Control 3	4.48	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
1430	165,000	1.9
358	55,500	2.2
89.4	15,200	2.6
22.3	4,060	2.4
5.59	1,100	1.5
1.40	333	2.8
0.349	130	5.0
0.00	64	5.7

## CERTIFICATE OF ANALYSIS

Page 2 of 2

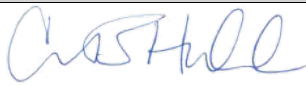
### Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-1 $\beta$ Plate	Z0047082	2-8°C	30 Sep 2021
Anti-ms IL-1 $\beta$ Antibody	D0081223	2-8°C	31 May 2022
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080168	2-8°C	31 May 2023
Diluent 41	M0410027	$\leq$ -10°C	31 Aug 2021
Diluent 45	M0450026	$\leq$ -10°C	30 Jun 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 (mouse) Control 1	Not Kit Specific
Proinflammatory Panel 1 (mouse) Control 2	Not Kit Specific
Proinflammatory Panel 1 (mouse) 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	Carissa Hubbard	Quality		19 JUN 2019