

## CERTIFICATE OF ANALYSIS

Page 1 of 2

**Product Description:** V-PLEX® and V-PLEX Plus NHP IL-1 $\beta$  Kit  
**Kit Catalog Numbers:** K156QP-Series  
**Kit Lot Number:** K0081141  
**Expiration Date:** 28 Feb 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	NHP 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	NHP IL-1 $\beta$	2.14	80% - 120%	< 20%	375	80% - 120%	< 20%

### Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	NHP IL-1 $\beta$	Calibrator	574	pg/mL
		Proinflammatory Panel 1 (human) Control 1	203	pg/mL
		Proinflammatory Panel 1 (human) Control 2	47.1	pg/mL
		Proinflammatory Panel 1 (human) Control 3	10.9	pg/mL

Controls tested may not be the same lot as shipped.

The calibrator and controls for this assay contain a human recombinant cytokine calibrator sequence that is highly homologous to the non-human primate cytokine. This assay has been verified to cross-react with homologs of rhesus and cynomolgus monkeys. See Product Insert for experimental data.

### Standard Curve Data:

NHP IL-1 $\beta$ - Spot 1		
Conc. pg/mL	ECL Signal	%CV
574	720,000	2.5
144	177,000	2.5
35.9	45,300	3.7
8.97	11,100	2.8
2.24	2,870	3.0
0.561	805	2.1
0.140	305	5.4
0.00	246	72.0

## CERTIFICATE OF ANALYSIS

Page 2 of 2

### Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-1 $\beta$ Plate	Z0046968	2-8°C	30 Apr 2021
Anti-hu IL-1 $\beta$ Antibody	D0081082	2-8°C	30 Sep 2021
Proinflammatory Panel 1 (human) Calibrator Blend	A0080180	2-8°C	30 Sep 2023
Diluent 2	M0020098	$\leq$ -10°C	28 Feb 2020
Diluent 3	M0030081	$\leq$ -10°C	30 Sep 2020
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
Review/Approval	Dawn Stollar	Quality Assurance		27 NOV 2018