

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human IL-1 β Kit

Kit Catalog Numbers: K151QP-series

Kit Lot Number: K0080942

Expiration Date: 31 Jul 2019

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	hu IL-1 β	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	hu IL-1 β	2.14	80% - 120%	< 20%	375	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu IL-1 β	Calibrator	566	pg/mL
		Proinflammatory Panel 1 (human) Control 1	215	pg/mL
		Proinflammatory Panel 1 (human) Control 2	52	pg/mL
		Proinflammatory Panel 1 (human) Control 3	11	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

hu IL-1 β - Spot 1		
Conc. pg/mL	ECL Signal	%CV
566	1,130,000	1.9
142	304,000	2.0
35.4	78,200	1.9
8.84	20,000	2.0
2.21	5,250	2.0
0.553	1,480	1.6
0.138	547	2.2
0.00	238	13.0

CERTIFICATE OF ANALYSIS

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-1 β Plate	Z0046726	2-8°C	30 Jun 2020
Anti-hu IL-1 β Antibody	D0080915	2-8°C	31 Jul 2019
Proinflammatory Panel 1 (human) Calibrator Blend	A0080147	2-8°C	31 Aug 2022
Diluent 2	M0020096	\leq -10°C	30 Nov 2019
Diluent 3	M0030077	\leq -10°C	30 Nov 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), 100 mL	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	<i>Dawn Stollar</i>	06 APR 2018