

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, K156QP-Series

Kit Lot Number: K0081511

Expiration Date: 31 Jan 2022

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 β	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 β	2.14	80% - 120%	< 20%	375	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	IL-1 β	Calibrator	610	pg/mL
		Proinflammatory Panel 1 (human) Control 1	235	pg/mL
		Proinflammatory Panel 1 (human) Control 2	55.9	pg/mL
		Proinflammatory Panel 1 (human) Control 3	11.9	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

IL-1 β - Spot 1		
Conc. pg/mL	ECL Signal	%CV
610	769,000	4.1
153	195,000	2.3
38.1	49,400	1.7
9.53	12,400	1.8
2.38	3,110	1.7
0.596	868	1.6
0.149	296	4.5
0.00	103	7.5

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-1 β Plate	Z0047356	2-8°C	31 Dec 2022
Anti-hu IL-1 β Antibody	D0081128	2-8°C	31 Jan 2022
Proinflammatory Panel 1 (human) Calibrator Blend	A0080242	2-8°C	31 Mar 2025
Diluent 2	M0020109	\leq -10°C	28 Feb 2022
Diluent 3	M0030096	\leq -10°C	30 Jun 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
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	Karen Hamilla	Quality		07 AUG 2020