

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human IL-1β Kit
Kit Catalog Numbers: K151QP-Series, K156QP-Series
Kit Lot Number: K0081527
Expiration Date: 31 Oct 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-1β	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

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

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	IL-1β	Calibrator	583	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-1β - Spot 1		
Conc. pg/ml	ECL Signal	%CV
583	736,000	2.6
146	186,000	1.6
36.4	46,700	1.6
9.11	11,600	2.0
2.28	2,900	1.9
0.569	804	2.6
0.142	289	3.1
0.00	122	7.7

CERTIFICATE OF ANALYSIS


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-1 β Plate	Z0047371	2-8°C	31 Jan 2023
Anti-hu IL-1 β Antibody	D0081380	2-8°C	31 Mar 2023
Proinflammatory Panel 1 (human) Calibrator blend	A0080262	2-8°C	30 Jun 2025
Diluent 2	M0020113	\leq -10°C	30 Nov 2022
Diluent 3	M0030100	\leq -10°C	31 Oct 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), 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	Karen Hamilla	Quality		22 SEP 2020