

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Mouse IL-1 β Kit
Kit Catalog Numbers: K152QP-Series
Kit Lot Number: K0081538
Expiration Date: 31 Dec 2022

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 β	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 β	0.720	75% - 125%	< 25%	1030	75% - 125%	< 25%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	IL-1 β	Calibrator	1530	pg/mL
		Proinflammatory Panel 1 (mouse) Control 1	722	pg/mL
		Proinflammatory Panel 1 (mouse) Control 2	52	pg/mL
		Proinflammatory Panel 1 (mouse) Control 3	4.32	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
1530	253,000	1.6
383	74,700	2.1
95.6	19,700	1.1
23.9	4,950	2.8
5.98	1,290	1.5
1.49	368	2.2
0.374	148	3.7
0.00	73	6.4

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-1 β Plate	Z0047381	2-8°C	28 Feb 2023
Anti-ms IL-1 β Antibody	D0081384	2-8°C	31 May 2023
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080252	2-8°C	30 Apr 2025
Diluent 41	M0410034	\leq -10°C	31 Dec 2022
Diluent 45	M0450035	\leq -10°C	31 Jan 2023
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	Karen Hamilla	Quality		06 OCT 2020