

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: K0081351
Expiration Date: 28 Feb 2021

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	1670	pg/mL
		Proinflammatory Panel 1 (mouse) Control 1	741	pg/mL
		Proinflammatory Panel 1 (mouse) Control 2	53.8	pg/mL
		Proinflammatory Panel 1 (mouse) Control 3	4.48	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
1670	376,000	2.3
418	107,000	2.1
104	28,000	2.2
26.1	7,170	1.8
6.52	1,830	1.5
1.63	501	2.1
0.408	168	2.2
0.00	63	7.8

CERTIFICATE OF ANALYSIS

Components:


Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-1β Plate	Z0047204	2-8°C	31 Mar 2022
Anti-ms IL-1B Antibody	D0081336	2-8°C	31 Jan 2023
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080075	2-8°C	28 Feb 2021
Diluent 41	M0410029	≤ -10°C	31 Dec 2021
Diluent 45	M0450030	≤ -10°C	28 Feb 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 (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		03 DEC 2019