

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: K0080973
Expiration Date: 31 May 2019

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	ms 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	ms IL-1 β	0.720	75% - 125%	< 25%	1030	75% - 125%	< 25%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	ms IL-1 β	Calibrator	1500	pg/mL
		Proinflammatory Panel 1 (mouse) Control 1	663	pg/mL
		Proinflammatory Panel 1 (mouse) Control 2	47.9	pg/mL
		Proinflammatory Panel 1 (mouse) Control 3	4.18	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

ms IL-1 β - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1500	247,000	2.1
375	73,600	3.1
93.8	20,700	4.4
23.4	5,240	3.2
5.86	1,430	4.5
1.46	397	4.5
0.366	152	3.1
0.00	69	6.2

CERTIFICATE OF ANALYSIS

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-1 β Plate	Z0046780	2-8°C	31 Aug 2020
Anti-ms IL-1 β Antibody	D0080852	2-8°C	31 May 2019
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080043	2-8°C	31 Jul 2019
Diluent 41	M0410022	\leq -10°C	31 Jul 2020
Diluent 45	M0450020	\leq -10°C	31 May 2020
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	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	17 APR 2018