

CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus Mouse IL-1β Kit
Kit Catalog Numbers: K152QP-series
Kit Lot Number: K0080730
Expiration Date: 28 Feb 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-1β	80% - 120%	< 20%	80% - 130%	< 25%	80% - 130%	< 25%	80% - 130%	< 25%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	ms IL-1β	0.720	75% - 125%	< 25%	1030	75% - 125%	< 25%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	ms IL-1β	Calibrator	1460	pg/mL
		Proinflammatory Panel 1 (mouse) Control 1	623	pg/mL
		Proinflammatory Panel 1 (mouse) Control 2	50.3	pg/mL
		Proinflammatory Panel 1 (mouse) Control 3	4.41	pg/mL

Controls tested may not be the same lot as shipped.

CERTIFICATE OF ANALYSIS

Page 2 of 2

Standard Curve Data:

ms IL-1 β - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1460	202,000	1.7
365	60,100	2.6
91.3	16,900	3.9
22.8	4,540	2.0
5.70	1,230	2.2
1.43	353	1.6
0.356	139	3.7
0.00	61	3.6


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IL-1 β Plate	Z0046405	2-8°C	31 Aug 2019
Mouse IL-1 β Antibody	D0080852	2-8°C	31 May 2019
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080060	2-8°C	31 Jul 2020
Diluent 41	M0410015	\leq -10°C	30 Jun 2019
Diluent 45	M0450012	\leq -10°C	28 Feb 2019
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		29 MAR 2017