

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Mouse IFN- γ Kit
Kit Catalog Numbers: K152QO-Series
Kit Lot Number: K0080792
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 IFN- γ	75% - 125%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	ms IFN- γ	0.390	75% - 125%	< 25%	570	75% - 125%	< 25%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	ms IFN- γ	Calibrator	933	pg/mL
		Proinflammatory Panel 1 (Mouse) Control 1	280	pg/mL
		Proinflammatory Panel 1 (Mouse) Control 2	21.5	pg/mL
		Proinflammatory Panel 1 (Mouse) Control 3	1.88	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

ms IFN- γ - Spot 1		
Conc. pg/mL	ECL Signal	%CV
933	392,000	1.9
233	92,200	4.7
58.3	22,500	2.7
14.6	5,590	1.8
3.64	1,440	2.4
0.911	410	5.1
0.228	155	2.2
0.00	68	5.2

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse IFN- γ Plate	Z0046483	2-8°C	30 Sep 2019
Anti-ms IFN- γ Antibody	D0080851	2-8°C	31 May 2019
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080110	2-8°C	31 Dec 2021
Diluent 41	M0410016	$\leq -10^{\circ}\text{C}$	31 Aug 2019
Diluent 45	M0450016	$\leq -10^{\circ}\text{C}$	30 Sep 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	<i>Dawn Stollar</i>	15 MAY 2017