

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-Plex and V- Plex® Plus Mouse KC/GRO Kit
Kit Catalog Numbers: K152QT-Series
Kit Lot Number: K0081098
Expiration Date: 30 Sep 2020

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 KC/GRO	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	ms KC/GRO	3.29	75% - 125%	< 25%	1230	75% - 125%	< 25%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	ms KC/GRO	Calibrator	1870	pg/mL
		Proinflammatory Panel 1 (mouse) Control 1	695	pg/mL
		Proinflammatory Panel 1 (mouse) Control 2	116	pg/mL
		Proinflammatory Panel 1 (mouse) Control 3	17.7	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

ms KC/GRO - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1870	802,000	1.9
468	176,000	2.4
117	37,400	2.0
29.2	9,040	1.1
7.30	2,120	1.0
1.83	599	2.2
0.457	205	2.9
0.00	74	5.4

CERTIFICATE OF ANALYSIS


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Mouse KC/GRO Plate	Z0046924	2-8°C	31 Mar 2021
Anti-ms KC/GRO Antibody	D0081152	2-8°C	28 Feb 2022
Proinflammatory Panel 1 (mouse) Calibrator Blend	A0080152	2-8°C	30 Nov 2022
Diluent 41	M0410024	≤ -10°C	30 Sep 2020
Diluent 45	M0450021	≤ -10°C	30 Sep 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		28 NOV 2018