

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Rat IL-6 Kit

Kit Catalog Numbers: K153QX-Series

Kit Lot Number: K0080961

Expiration Date: 30 Sep 2019

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Proinflammatory Panel 1 (rat) Control 1		Proinflammatory Panel 1 (rat) Control 2		Proinflammatory Panel 1 (rat) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	rt IL-6	80% - 120%	< 20%	75% - 125%	< 25%	75% - 125%	< 25%	75% - 125%	< 25%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	rt IL-6	96.9	75% - 125%	< 25%	8550	75% - 125%	< 25%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	rt IL-6	Calibrator	13300	pg/mL
		Proinflammatory Panel 1 (rat) Control 1	10600	pg/mL
		Proinflammatory Panel 1 (rat) Control 2	3480	pg/mL
		Proinflammatory Panel 1 (rat) Control 3	1130	pg/mL

Controls tested may not be the same lot as shipped.

Standard Curve Data:

rt IL-6 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
13300	85,200	3.9
3330	17,700	4.7
831	4,210	3.8
208	1,140	3.6
52.0	381	2.2
13.0	185	3.1
3.25	140	3.9
0.00	135	19.0

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Rat IL-6 Plate	Z0046763	2-8°C	31 Jul 2020
Anti-rat IL-6 Antibody	D0081029	2-8°C	29 Feb 2020
Proinflammatory Panel 1 (rat) Calibrator Blend	A0080100	2-8°C	30 Sep 2021
Blocker H	R0100015	Room Temperature	30 Sep 2019
Diluent 42	M0420010	≤ -10°C	31 Jul 2020
Diluent 40	M0400011	≤ -10°C	31 Aug 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), 1000mL	Not Kit Specific
Proinflammatory Panel 1 (rat) Control 1	Not Kit Specific
Proinflammatory Panel 1 (rat) Control 2	Not Kit Specific
Proinflammatory Panel 1 (rat) 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		02 JUL 2018