

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human ICAM-1 Kit
Kit Catalog Numbers: K151SUD-series
Kit Lot Number: K0080926
Expiration Date: 30 Sep 2019

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Vascular Injury Control 1		Vascular Injury Control 2		Vascular Injury Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	hu ICAM-1	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	hu ICAM-1	6.40	80% - 120%	< 20%	32700	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu ICAM-1	20X Calibrator	968000	pg/mL
		Vascular Injury Control 1	16300	pg/mL
		Vascular Injury Control 2	1720	pg/mL
		Vascular Injury Control 3	209	pg/mL

Controls tested may not be the same lot as shipped.

Concentrations provided in this table refer to the stock concentration of the calibrators. For top of the standard curve concentrations, please refer to the Standard Curve Data

Standard Curve Data:

hu ICAM-1 - Spot 1		
Conc. pg/mL	ECL Signal	%CV
48400	664,000	6.0
9680	147,000	4.5
1940	29,100	6.6
387	5,770	4.7
77.4	1,110	5.5
15.5	242	4.1
3.10	75	8.2
0.00	39	12.0

Components:


CERTIFICATE OF ANALYSIS

Description	Lot Number	Storage Temperature	Expiration Date
Human ICAM-1 Plate	Z0056115	2-8°C	31 May 2020
Anti-hu ICAM-1 Antibody	D0080981	2-8°C	30 Sep 2019
Vascular Injury Panel 2 (human) Calibrator Blend	A0080132	≤ -70°C	30 Sep 2019
Diluent 101	M1010029	≤ -10°C	31 Mar 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
Vascular Injury Control 1	Not Kit Specific
Vascular Injury Control 2	Not Kit Specific
Vascular Injury 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		15 DEC 2017