

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Human CRP Kit

Kit Catalog Numbers: K151ST Series

Kit Lot Number: K0080900

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 CRP	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	hu CRP	27.6	80% - 120%	< 20%	49600	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	hu CRP	20X Calibrator	4040000	pg/mL
		Vascular Injury Control 1	24300	pg/mL
		Vascular Injury Control 2	4900	pg/mL
		Vascular Injury Control 3	604	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 CRP - Spot 1		
Conc. pg/mL	ECL Signal	%CV
202000	963,000	3.1
40400	364,000	3.7
8080	84,100	3.8
1620	17,200	2.6
323	3,510	3.6
64.6	725	4.3
12.9	164	4.7
0.00	25	22.0

Components:

CERTIFICATE OF ANALYSIS

Page 2 of 2

Description	Lot Number	Storage Temperature	Expiration Date
Human CRP Plate	Z0056078	2-8°C	31 Mar 2020
Anti-hu CRP Antibody	D0081021	2-8°C	31 Jan 2020
Vascular Panel 2(Human) Calibrator Blend	A0080132	≤ -70°C	30 Sep 2019
Diluent 101	M1010028	≤ -10°C	29 Feb 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
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	<i>Dawn Stollar</i>	10 NOV 2017