

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: K0081152
Expiration Date: 31 Mar 2021

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	3680000	pg/mL
		Vascular Injury Control 1	25100	pg/mL
		Vascular Injury Control 2	5060	pg/mL
		Vascular Injury Control 3	639	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
184000	858,000	1.6
36800	310,000	2.5
7360	67,500	3.9
1470	13,200	2.4
294	2,530	2.7
58.9	529	2.9
11.8	130	6.1
0.00	24	15.0

CERTIFICATE OF ANALYSIS

Page 2 of 2


Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human CRP Plate	Z0056409	2-8°C	30 Apr 2021
Ant-hu CRP antibody	D0081221	2-8°C	31 May 2022
Vascular Injury Panel 2 (human) Calibrator Blend	A0080184	≤ -70°C	31 Mar 2021
Diluent 101	M1010033	≤ -10°C	30 Apr 2021
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
MSD Wash Buffer (20x), 100mL	Not Kit Specific
Plate Seals	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	Karen Hamilla	Quality		26 SEP 2019