

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus Vascular Injury Panel 2 (human) Kit
Kit Lot Number: K0081551
Expiration Date: 31 Jul 2021
Kit Catalog Numbers: K15198-Series; K151SS-Series, K151SR-Series; K151SU-Series

Precision and Accuracy Specifications:

		Calibrator		Vascular Injury Control 1		Vascular Injury Control 2		Vascular Injury Control 3	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	SAA	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
2	CRP	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
3	VCAM-1	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%
4	ICAM-1	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	SAA	54.0	80% - 120%	< 20%	138000	80% - 120%	< 20%
2	CRP	27.6	80% - 120%	< 20%	49600	80% - 120%	< 20%
3	VCAM-1	37.6	80% - 120%	< 20%	32000	80% - 120%	< 20%
4	ICAM-1	15.0	80% - 120%	< 20%	32700	80% - 120%	< 20%

Non-Specific Binding Specifications:

Spot	Name	NSB Spec
1	SAA	< 1.0%
2	CRP	< 1.0%
3	VCAM-1	< 1.0%
4	ICAM-1	< 1.0%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	SAA	4940000	pg/mL
2	CRP	3760000	pg/mL
3	VCAM-1	994000	pg/mL
4	ICAM-1	998000	pg/mL

CERTIFICATE OF ANALYSIS

Page 2 of 2

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Vascular Injury Panel 2 (human) Plate	Z0056794	2-8°C	28 Feb 2023
Anti-hu SAA Antibody	D0081403	2-8°C	31 Jul 2023
Anti-hu CRP Antibody	D0081563	2-8°C	31 Mar 2024
Anti-hu VCAM-1 Antibody	D0081564	2-8°C	31 Mar 2024
Anti-hu ICAM-1 Antibody	D0081565	2-8°C	31 Mar 2024
Vascular Injury Panel 2 (human) Calibrator Blend	A0080249	≤ -70°C	30 Nov 2022
Diluent 101	M1010030	≤ -10°C	31 Jul 2021

Additional components may be provided to execute the assay. Additional components are not lot-specific.

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		08 JAN 2021