



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX® and V-PLEX Plus Human VEGF Kit
Lot Number: K0082958
Expiration Date: 30 SEP 2027
Catalog Number: K151RH-Series, K156RH-Series

Lot-Specific Components:

The kit expiration date provided in the COA is determined by the component with the earliest expiration date. As the kit is supplied with different catalog products, not all configurations include all the components. In such cases, the applicable expiration date is determined by the shortest expiry among the components included in the kit that is received.

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu VEGF Antibody	D0082352	2-8°C	30 APR 2028
Cytokine Panel 1 (human) Calibrator Blend	A0080454	2-8°C	31 DEC 2030
Human VEGF-A 96-Well Small Spot SECTOR Plate	Z0057706	2-8°C	30 NOV 2028
Diluent 3	M0030148	≤ -10°C	31 JUL 2028
Diluent 43	M0430059	≤ -10°C	30 SEP 2027

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

Precision, Accuracy, and Sensitivity Specifications:

		Calibrator		Cytokine Panel 1 (human) Control 1		Cytokine Panel 1 (human) Control 2		Cytokine Panel 1 (human) Control 3	
Spot	Name	Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	VEGF-A	80% - 120%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%

		LLOQ			ULOQ		
Spot	Name	Value	Recovery	CV	Value	Recovery	CV
1	VEGF-A	7.70	80% - 120%	< 20%	562	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	VEGF-A	1150	pg/mL

Dilution to top of Curve: 1X

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		16 JUL 2026