



Meso Scale Diagnostics, LLC.

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human IL-8 Kit
Kit Lot Number: K0081962
Expiration Date: 30 Sep 2023
Kit Catalog Numbers: K151RA-Series; K156RA-Series

Precision and Accuracy Specifications:

Spot	Name	Calibrator		Proinflammatory Panel 1 (human) Control 1		Proinflammatory Panel 1 (human) Control 2		Proinflammatory Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-8	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-8	0.591	80% - 120%	< 20%	375	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	IL-8	601	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human IL-8 Plate	Z0047661	2-8°C	31 Jul 2024
Anti-hu IL-8 Antibody	D0081700	2-8°C	30 Nov 2024
Proinflammatory Panel 1 (human) Calibrator Blend	A0080299	2-8°C	30 Jun 2026
Diluent 2	M0020116	≤ -10°C	30 Sep 2023
Diluent 3	M0030111	≤ -10°C	31 May 2024

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		29 APR 2022