

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human MIP-1β Kit
Kit Catalog Numbers: K151NR-Series, K156NR-Series
Kit Lot Number: K0081588
Expiration Date: 30 Nov 2022

Precision and Accuracy Specifications:

Spot	Name	Calibrator		Chemokine Panel 1 (human) Control 1		Chemokine Panel 1 (human) Control 2		Chemokine Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	MIP-1β	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	MIP-1β	1.88	80% - 120%	< 20%	750	80% - 120%	< 20%

Stock Concentration of Calibrator:


Spot	Name	Conc.	Units
1	MIP-1β	1190	pg/mL

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human MIP-1β Plate	Z0047416	2-8°C	30 Apr 2023
Anti-hu MIP-1β Antibody	D0081312	2-8°C	31 Jan 2023
Chemokine Panel 1 (human) Calibrator Blend	A0080266	2-8°C	31 Aug 2025
Diluent 43	M0430040	≤ -10°C	30 Nov 2022
Diluent 3	M0030101	≤ -10°C	31 Dec 2022

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		15 DEC 2020