

CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: V-PLEX® and V-PLEX Plus Human MIP-1 β Kit
Kit Lot Number: K0081633
Expiration Date: 31 Jul 2022
Kit Catalog Numbers: K151NR-Series, K156NR-Series

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-1 β	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%	80% - 120%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	MIP-1 β	2.82	80% - 120%	< 20%	520	80% - 120%	< 20%

Stock Concentration of Calibrator:

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

Lot-Specific Components:


Description	Lot Number	Storage Temperature	Expiration Date
Human MIP-1 β Plate	Z0047455	2-8°C	30 Jun 2023
Anti-hu MIP-1 β Antibody	D0081181	2-8°C	31 Jul 2022
Chemokine Panel 1 (human) Calibrator Blend	A0080266	2-8°C	31 Aug 2025
Diluent 43	M0430037	\leq -10°C	31 Aug 2022
Diluent 3	M0030103	\leq -10°C	30 Apr 2023

Additional Comments:

Assay	Spot	Specification	Comment
MIP-1 β	1	LLOQ Conc.	Did not meet specifications at 1.88 pg/mL; LLOQ reported at 2.82 pg/mL.

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		23 JUN 2021