

CERTIFICATE OF ANALYSIS

Page 1 of 2

Product Description: V-PLEX® and V-PLEX Plus NHP MIP-1β Kit
Kit Catalog Numbers: K156NR-Series
Kit Lot Number: K0080944
Expiration Date: 30 Sep 2019

Acceptance Data/ Percent Recovery 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	NHP 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	NHP MIP-1β	2.27	80% - 120%	< 20%	750	80% - 120%	< 20%

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	NHP MIP-1β	Calibrator	1110	pg/mL
		Chemokine Panel 1 (human) Control 1	734	pg/mL
		Chemokine Panel 1 (human) Control 2	116	pg/mL
		Chemokine Panel 1 (human) Control 3	17.1	pg/mL

Controls tested may not be the same lot as shipped.

The calibrator and controls for this assay contain a human recombinant cytokine calibrator sequence that is highly homologous to the non-human primate cytokine. This assay has been verified to cross-react with homologs of rhesus and cynomolgus monkeys. See Product Insert for experimental data.

Standard Curve Data:

NHP MIP-1β - Spot 1		
Conc. pg/mL	ECL Signal	%CV
1110	511,000	2.8
278	114,000	1.9
69.4	21,100	2.2
17.3	3,740	2.9
4.34	925	3.4
1.08	266	3.6
0.271	113	3.6
0.00	63	6.3

CERTIFICATE OF ANALYSIS

Page 2 of 2

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Human MIP-1 β Plate	Z0046728	2-8°C	30 Jun 2020
Anti-hu MIP-1 β Antibody	D0080898	2-8°C	30 Sep 2019
Chemokine Panel 1 (human) Calibrator Blend	A0080122	2-8°C	28 Feb 2022
Diluent 43	M0430021	\leq -10°C	30 Sep 2019
Diluent 3	M0030076	\leq -10°C	30 Sep 2019
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Additional Components Included in V-PLEX Plus	
Plate Seals	Not Kit Specific
Wash Buffer (20X), 100mL	Not Kit Specific
Chemokine Panel 1 (human) Control 1	Not Kit Specific
Chemokine Panel 1 (human) Control 2	Not Kit Specific
Chemokine Panel 1 (human) Control 3	Not Kit Specific

Statement:

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

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	21 FEB 2018