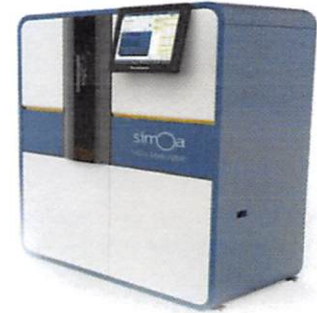


Product Number:	101160
Lot Number:	501741
Expiration:	13-Mar-2020
Platform(s):	HD-1 & SR-X



Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent †	101263	908701
Detector Reagent	101264	908702
SBG Reagent	101296	907212
Sample Diluent	103329	907211
Calibrators	101299	81720/826802
Controls	101302/101316	81721
SBG Diluent (SR-X)††	103158	N/A ¹
RGP Reagent	103159	N/A ¹

† Follow the Simoa™ Cytokine 3-Plex A Advantage Kit Instructions (KI-0011 - HD1/KI-0101 - SR-X) to prepare the Bead Reagent.
†† Specific for use on the SR-X platform - refer to the Simoa™ Cytokine 3-Plex A SR-X Kit Instructions (KI-0101).

¹ Reagents are not Kit Lot Specific

Release Materials

IL-6

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	81721	0.942	0.717-1.08
Control 2	81721	26.1	19.7-29.5
Panel 2	MF-062818-02	0.983	0.776-1.16
Panel 3	MF-062818-03	1.65	1.40-2.10
Panel 4	MF-062818-04	13.7	10.2-15.3

IL-10

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	81721	0.683	0.536-0.803
Control 2	81721	16.7	10.8-18.8
Panel 2	MF-062818-02	14.1	9.27-16.0
Panel 3	MF-062818-03	0.594	0.419-0.768
Panel 4	MF-062818-04	1.34	1.00-1.50

Ranges shown are generated internally for new lot release only.

Customer should generate their own control ranges.

TNFα

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	81721	1.98	1.52-2.28
Control 2	81721	51.9	39.0-58.5
Panel 2	MF-062818-02	2.06	1.54-2.31
Panel 3	MF-062818-03	38.2	31.6-47.4
Panel 4	MF-062818-04	6.78	5.40-8.10

Review/Approval

Patrizia Stadler

Assoc. Director, QC

P. Stadler 22 Apr 2019

Name

Title

Signature/Date