

Product Number:	101319
Lot Number:	501660
Expiration:	29-Jun-2019
Platform(s):	HD-1



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

Component	Part Number	Lot Number
Bead Reagent	101265	826815
Detector Reagent	101266	826814
SBG Reagent (HD-1)†	101297	823312
Sample Diluent	103330	826206
Calibrators	101300	80520
Controls	101303 & 101317	80521
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials

IL-6

	Lot	Result (pg/mL)	Historical Range (pg/mL)
Control 1	80521	1.05	0.838-1.26
Control 2	80521	29.4	22.9-34.4
Panel 2	MF-062818-02	0.927	0.782-1.17
Panel 3	MF-062818-03	1.55	1.27-1.91
Panel 4	MF-062818-04	13.1	10.8-16.2

IL-17A

	Lot	Result (pg/mL)	Historical Range (pg/mL)
Control 1	80521	0.803	0.651-0.976
Control 2	80521	17.4	13.6-20.4
Panel 1	MF-062818-02	17.1	14.0-21.1
Panel 3	MF-062818-03	1.52	1.18-1.78
Panel 4	MF-062818-04	2.07	1.67-2.51

TNFα

	Lot	Result (pg/mL)	Historical Range (pg/mL)
Control 1	80521	2.00	1.57-2.36
Control 2	80521	49.8	39.2-58.7
Panel 1	MF-062818-02	2.02	1.51-2.27
Panel 3	MF-062818-03	40.1	33.7-50.6
Panel 4	MF-062818-04	7.11	5.63-8.44

Ranges shown are generated internally for new lot release only.

Customer should generate their own target ranges.

Review/Approval	Patrizia Stadler	QC Assoc. Director	<i>P. Stadler</i>	15 Feb 2019
	Name	Title	Signature/Date	