

## Certificate of Analysis

Kit Name:	Simoa™ IL-17A 2.0 Kit
Product Number:	101599
Lot Number:	500895
Expiration:	1-Jun-2018

113 Hartwell Ave.  
Lexington MA.  
02421

[www.quanterix.com](http://www.quanterix.com)

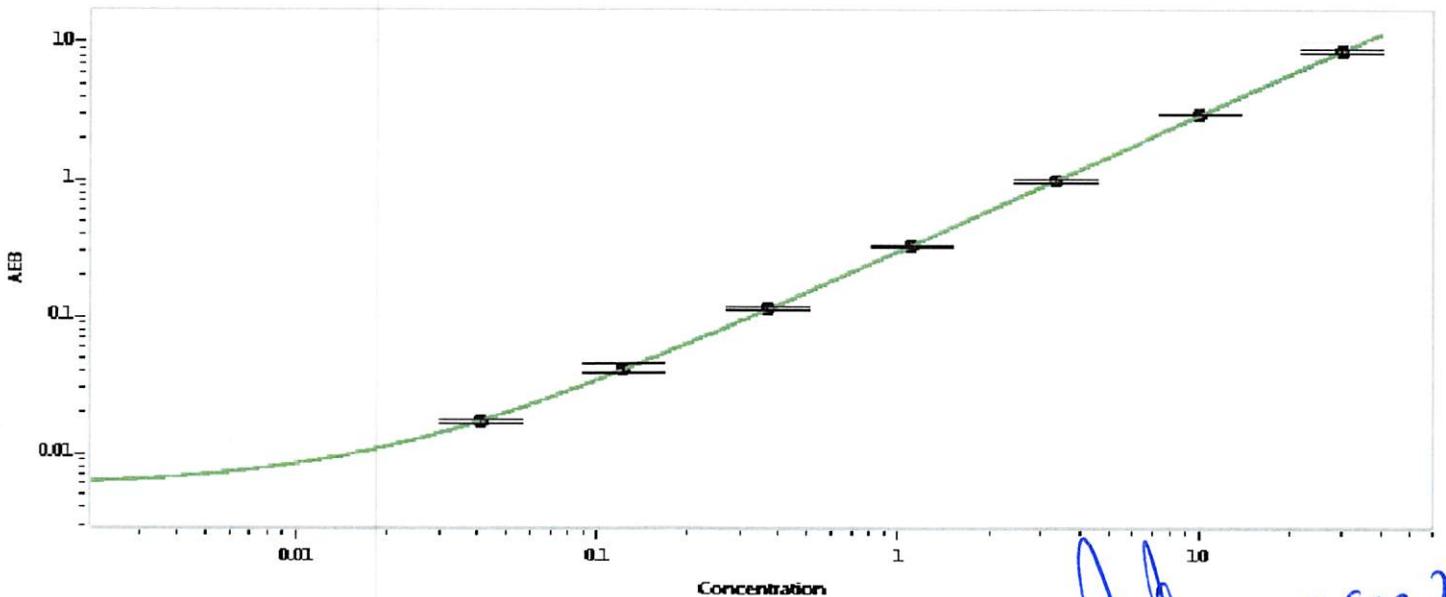
Component	Part Number	Lot Number
IL-17A 2.0 Bead Reagent	101601	72492
IL-17A 2.0 Detector Reagent	102012	72493
IL-17A 2.0 SBG Reagent	101602	72494
IL-17A 2.0 Sample Diluent	101603	72441
IL-17A 2.0 Calibrator Diluent	102014	72443
IL-17A Calibrator Concentrate	100778	70764
RGP Reagent	100030	1

<sup>1</sup> RGP is not Kit Lot Specific

Calibrator Concentrate	Assigned Value (pg/mL)
	636

Controls	Lot #s	Range (pg/mL)	Value at Release (pg/mL)
Control 1	52442	0.485-0.727	0.633
Control 2	52442	19.5-34.7	29.3

Patient Panels	Lot Number	Range (pg/mL)	Result (pg/mL)
1	MF-071615-01	0.325-0.513	0.440
2	MF-071615-02	5.58-8.37	8.14
3	MF-071615-03	1.68-2.52	2.38



James Drummey

Senior Director  
Manufacturing & Operations

*19-Sep-2017*

*[Handwritten Signature]*

### Review/Approval

Name

Title

Signature/Date

Statement: The above product was tested using Quanterix procedures and conforms to Quanterix's specifications. It is intended for Research Use Only. Not for use in Diagnostic Procedures.