

Certificate of Analysis

Kit Name:	Simoa™ Homebrew Assay Development Kit
Product Number:	101354
Lot Number:	500916
Expiration:	16-Aug-2018


113 Hartwell Ave.
Lexington MA.
02421
www.quanterix.com

Component	Part Number	Lot Number
Bead Diluent Buffer	100458	MF-081617-03
Bead Wash Buffer	101355	MF-081817-11
Bead Blocking Buffer	101356	MF-081617-09
Bead Conjugation Buffer	101357	MF-081617-06
Biotinylation Reaction Buffer	101358	MF-081817-06
Detector/Sample Diluent	101359	MF-081817-09
Beads	101360	MF-091517-07

Bead Count (beads/mL)	
Assigned Value	2.76X10 ⁹

James Drummey

Senior Director
Manufacturing & Operations

 22-Sep-2017

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.