

Kit Name:	Simoa Qualification Test
Product Number:	101537
Lot Number:	500968
Expiration:	26-Apr-2018

Component	Part Number	Lot Number
SQT Bead	101535	73041
SQT Detector	101533	73000
SQT SBG	101534	72291
SQT Controls	101543-101544	73060
RGP	100030	1

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

Bead Count	1.59X10 <sup>7</sup>
Bead % Monomeric	91%

Review/Approval

Sivasankari Murugan

QC Supervisor

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.