

|                        |             |
|------------------------|-------------|
| <b>Product Number:</b> | 101361      |
| <b>Lot Number:</b>     | 501415      |
| <b>Expiration:</b>     | 10-Jul-2019 |

| Component           | Part Number | Lot Number       | Concentration |
|---------------------|-------------|------------------|---------------|
| SBG Dilution Buffer | 100376      | 821410           | N/A           |
| SBG Concentrate     | 100439      | 819110           | 151 nM        |
| RGP Reagent         | 103159      | N/A <sup>1</sup> | N/A           |

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

**Review/Approval**

Sivasankari Murugan  
**Name**

QC Supervisor  
**Title**

 18 SEP 2018  
**Signature/Date**

**Statement:** The above product was tested using Quanterix procedures and conforms to Quanterix's specifications.