

<b>Product Number:</b>	101361
<b>Lot Number:</b>	501534
<b>Expiration:</b>	22-Aug-2019

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	829915	N/A
SBG Concentrate	100439	823409	155 nM
RGP Reagent	103159	N/A <sup>1</sup>	N/A

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

**Review/Approval**

Sivasankari Murugan

QC Supervisor

Name

Title

*Sivasankari Murugan* 06 DEC 2018

Signature/Date

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