

<b>Product Number:</b>	101361
<b>Lot Number:</b>	501445
<b>Expiration:</b>	10-Jul-2019

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	823703	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      QC Supervisor            05 OCT 2018  
Name      Title      Signature/Date

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