

<b>Product Number:</b>	101361
<b>Lot Number:</b>	501731
<b>Expiration:</b>	11-Mar-2020

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	907014	N/A
SBG Concentrate	103397	907013	50 nM
RGP Reagent	103159	N/A <sup>1</sup>	N/A

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

Review/Approval

Patrizia Stadler

Assoc. Director, QC

25 Mar 2019 P. Stadler

Name

Title

Signature/Date

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