

<b>Product Number:</b>	101361
<b>Lot Number:</b>	501679
<b>Expiration:</b>	19-Nov-2019

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	834903	N/A
SBG Concentrate	103397	832317	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

*P. Stadler* 25 Feb 2019

Name

Title

Signature/Date

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