

<b>Product Number:</b>	101361
<b>Lot Number:</b>	501589
<b>Expiration:</b>	29-Aug-2019

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	834903	N/A
SBG Concentrate	100439	824115	155 nM
RGP Reagent	103159	N/A <sup>1</sup>	N/A

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

Review/Approval	Patrizia Stadler Associate Director, QC	<i>P. Stadler</i> 24 JAN 2019
	Name	Title
		Signature/Date

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