

<b>Product Number:</b>	101361
<b>Lot Number:</b>	502442
<b>Expiration:</b>	2-May-2021

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	012309	N/A
SBG Concentrate	103397	013908	50 nM
RGP Reagent	103159	N/A <sup>1</sup>	N/A

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

Review/Approval

Shivani Goel  
Name

QC Manager  
Title

*Shivani Goel 29 Jun 2020*  
Signature/Date