

<b>Product Number:</b>	101361
<b>Lot Number:</b>	503230
<b>Expiration:</b>	28-Jun-2022

Component	Part Number	Lot Number	Concentration
<b>SBG Dilution Buffer</b>	100376	121714	N/A
<b>SBG Concentrate</b>	103397	120031	50 nM
<b>RGP Reagent</b>	103159	N/A <sup>1</sup>	N/A

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

Review/Approval

*Shivani Goel*  
Name

QC

Title

*Shivani Goel 30 Dec 2021*  
Signature/Date