

<b>Product Number:</b>	101361
<b>Lot Number:</b>	502797
<b>Expiration:</b>	15-Jan-2022

Component	Part Number	Lot Number	Concentration
<b>SBG Dilution Buffer</b>	100376	104728	N/A
<b>SBG Concentrate</b>	103397	101913	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

*Junpil Park*  
Name

QC Supervisor  
Title

*[Signature]* 11 Mar 2021  
Signature/Date