

<b>Product Number:</b>	101361
<b>Lot Number:</b>	502986
<b>Expiration:</b>	15-Mar-2022

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	112808	N/A
SBG Concentrate	103397	107425	50 nM
RGP Reagent	103159	N/A <sup>1</sup>	N/A

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

Review/Approval

Jungil Park  
Name

QC supervisor  
Title

[Signature] (11 June 2021)  
Signature/Date