

<b>Product Number:</b>	101361
<b>Lot Number:</b>	503058
<b>Expiration:</b>	1-Jun-2022

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	118214	N/A
SBG Concentrate	103397	116813	50 nM
RGP Reagent	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]* 16 Jul 2021

Signature/Date