

<b>Product Number:</b>	101361
<b>Lot Number:</b>	502099
<b>Expiration:</b>	22-Apr-2020

Component	Part Number	Lot Number	Concentration
<b>SBG Dilution Buffer</b>	100376	921119	N/A
<b>SBG Concentrate</b>	103397	911210	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

Patrizia Stadler

Assoc. Director, QC

*P. Stadler* 01 Nov 2019

Name

Title

Signature/Date