

<b>Product Number:</b>	101361
<b>Lot Number:</b>	502264
<b>Expiration:</b>	15-Jan-2021

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	001516	N/A
SBG Concentrate	103397	002714	50 nM
RGP Reagent	103159	N/A <sup>1</sup>	N/A

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

Review/Approval

Patrizia Stadler  
Name

Assoc. Director, QC  
Title

*p.stadler* *9Mar2020*  
Signature/Date