

Product Number:	101361
Lot Number:	503669
Expiration:	22-Feb-2023

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	217316	N/A
SBG Concentrate	103397	209718	50 nM
RGP Reagent	103159	N/A <sup>1</sup>	N/A

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

Review/Approval

*Sonia Khillan*  
Name

QC  
Title

*Sonia Khillan / 08 AUG 2022*  
Signature/Date