

Product Number:	101361
Lot Number:	503517
Expiration:	16-Feb-2022

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	208911	N/A
SBG Concentrate	103397	209717	50 nM
RGP Reagent	103159	N/A ¹	N/A

¹ RGP is not Kit Lot Specific

Review/Approval

Matt Smart
Name

QC
Title

[Signature] 09 Jun 2022
Signature/Date