

Product Number:	101361
Lot Number:	503691
Expiration:	23-May-2023

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	222007	N/A
SBG Concentrate	103397	215710	50 nM
RGP Reagent	103159	N/A ¹	N/A

¹ RGP is not Kit Lot Specific

Review/Approval

Sonia Khillan

Name

QC

Title

Sonia Khillan / 08SEP2022

Signature/Date