

Product Number:	101361
Lot Number:	503803
Expiration:	29-Aug-2023

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	302512	N/A
SBG Concentrate	103397	230622	50 nM
RGP Reagent	103159	N/A ¹	N/A

¹ RGP is not Kit Lot Specific

Review/Approval

Sonia Khillan

Name

QC Manager

Title

Sonia Khillan 08 FEB 2023

Signature/Date