

Product Number:	101361
Lot Number:	503968
Expiration:	14-Jun-2024

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	316508	N/A
SBG Concentrate	103397	315712	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 07SEP2023
Signature/Date