

Product Number:	101361
Lot Number:	503340
Expiration:	30-Aug-2022

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	133217	N/A
SBG Concentrate	103397	124215	50 nM
RGP Reagent	103159	N/A ¹	N/A

¹ RGP is not Kit Lot Specific

Review/Approval

Shivani Goel
Name

QC

Title

Shivani Goel 02 Feb 2022
Signature/Date