

Product Number:	101361
Lot Number:	502716
Expiration:	15-Dec-2021

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	100805	N/A
SBG Concentrate	103397	101116	50 nM
RGP Reagent	103159	N/A ¹	N/A

¹ RGP is not Kit Lot Specific

Review/Approval

Shivani Goel
Name

Manager, QC
Title

Shivani Goel 10 Feb 2021
Signature/Date