

Product Number:	101361
Lot Number:	502032
Expiration:	30-Jul-2020

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	921117	N/A
SBG Concentrate	103397	921713	50 nM
RGP Reagent	103159	N/A ¹	N/A

¹ RGP is not Kit Lot Specific

Review/Approval

Patrizia Stadler

Assoc. Director, QC

P. Stadler

19 Sep 2019

Name

Title

Signature/Date