

Product Number:	101361
Lot Number:	502322
Expiration:	27-Jan-2021

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	005225	N/A
SBG Concentrate	103397	002715	50 nM
RGP Reagent	103159	N/A ¹	N/A

¹ RGP is not Kit Lot Specific

Review/Approval

Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i>	14 Apr 2020
Name	Title	Signature	Date