

Product Number:	101361
Lot Number:	501850
Expiration:	15-Apr-2020

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	916108	N/A
SBG Concentrate	103397	910511	50 nM
RGP Reagent	103159	N/A ¹	N/A

¹ RGP is not Kit Lot Specific

Review/Approval

Patrizia Stadler
Name

Assoc. Director, QC
Title

P. Stadler 21 JUN 2019
Signature/Date