

Product Number:	101361
Lot Number:	501664
Expiration:	19-Nov-2019

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	834903	N/A
SBG Concentrate	103397	832316	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 14 Feb 2019
Signature/Date

Statement: The above product was tested using Quanterix procedures and conforms to Quanterix's specifications.