

Product Number:	101361
Lot Number:	501622
Expiration:	29-Aug-2019

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	834903	N/A
SBG Concentrate	100439	824116	155 nM
RGP Reagent	103159	N/A ¹	N/A

¹ RGP is not Kit Lot Specific

Review/Approval

Patrizia Stadler Assoc. Director, QC
Name Title

p. Stadler 30 Jan 2019
Signature/Date

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