

Product Number:	101361
Lot Number:	501845
Expiration:	11-Mar-2020

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	907014	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 Assoc. Director, QC
Name Title

P. Stadler 21 May 2019
Signature/Date

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