

Product Number:	101361
Lot Number:	501528
Expiration:	22-Aug-2019

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	829912	N/A
SBG Concentrate	100439	823409	155 nM
RGP Reagent	103159	N/A ¹	N/A

¹ RGP is not Kit Lot Specific

Review/Approval

Sivasankari Murugan

QC Supervisor

Name

Title

 06 DEC 2018

Signature/Date

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