

Product Number:	101361
Lot Number:	501217
Expiration:	3-Oct-2018

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	MF-013018-15	
SBG Concentrate	100439	72765	165 nM
RGP Reagent	103159	N/A ¹	

¹ RGP is not Kit Lot Specific

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title


Signature/Date

16 APR 2018

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