

Product Number:	101361
Lot Number:	501156
Expiration:	19-Dec-2018

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	MF-121917-13	
SBG Concentrate	100439	80373	119 nM
RGP Reagent	103159	N/A ¹	

¹ RGP is not Kit Lot Specific

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title


Signature/Date 27 FEB 2018

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