


Product Number:	101361
Lot Number:	501466
Expiration:	6-Aug-2019

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	823704	N/A
SBG Concentrate	100439	821805	144 nM
RGP Reagent	103159	N/A ¹	N/A

¹ RGP is not Kit Lot Specific

Review/Approval	Sivasankari Murugan	QC Supervisor	
	Name	Title	Signature/Date

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