


Product Number:	101361
Lot Number:	501512
Expiration:	10-Jul-2019

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	823704	N/A
SBG Concentrate	100439	819110	151 nM
RGP Reagent	103159	N/A ¹	N/A

¹ RGP is not Kit Lot Specific

	James Drummey	
	Sr. Director, Manufacturing & Operations	
	on behalf of:	
Review/Approval	Sivasankari Murugan	QC Supervisor
	Name	Title
		
		26-Nov-2018
		Signature/Date

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