


Product Number:	101361
Lot Number:	501248
Expiration:	4-Apr-2019

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	MF-041518-05	
SBG Concentrate	100439	80946	164 nM
RGP Reagent	103159	N/A ¹	

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.