

Product Number:	101361
Lot Number:	501128
Expiration:	28-Nov-2018

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	MF-121917-13	
SBG Concentrate	100439	73328	128 nM
RGP Reagent	103159	N/A <sup>1</sup>	

<sup>1</sup> RGP is not Kit Lot Specific

Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

  
Signature/Date

14 FEB 2018

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