

Product Number:	101361
Lot Number:	501235
Expiration:	2-Apr-2019
Storage:	2-8°C

Component	Part Number	Lot Number	Concentration
SBG Dilution Buffer	100376	MF-041518-05	
SBG Concentrate	100439	80925	144 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

 30 APR 2018  
Signature/Date

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