

Product Number:	101361
Lot Number:	501331
Expiration:	23-May-2019

Component	Part Number	Lot Number
SBG Dilution Buffer	100376	MF-052318-12
SBG Concentrate	100439	815603
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

 19/05/2018  
Signature/Date

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