

Product Number:	101735
Lot Number:	501334
Expiration:	23-May-2019

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

SBG Concentrate Concentration (nM)	159
------------------------------------	-----

Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

  
Signature/Date

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