

Product Number:	101735
Lot Number:	501547
Expiration:	29-Aug-2019

Component	Item Number	Lot Number
SBG Concentrate	100439	824115
SBG Dilution Buffer	100376	829916

SBG Concentrate Concentration (nM)	155
---------------------------------------	-----

Review/Approval

Sivasankari Murugar
Name

QC Supervisor
Title

 12DEC2018
Signature/Date

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