

Product Number:	101735
Lot Number:	501467
Expiration:	6-Aug-2019

Component	Item Number	Lot Number
SBG Concentrate	100439	821805
SBG Dilution Buffer	100376	823704

SBG Concentrate Concentration (nM)	144
------------------------------------	-----

Review/Approval

Sivasankari Murugar QC Supervisor
Name Title

 23 OCT 2018
Signature/Date

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