

Product Number:	102295
Lot Number:	501395
Expiration:	10-Jul-2019

Component	Item Number	Lot Number
SBG Concentrate	100439	819109

SBG Concentrate Concentration (nM)	151
---------------------------------------	-----

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

 05 SEP 2018
Signature/Date

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