

Product Number:	102295
Lot Number:	501505
Expiration:	6-Aug-2019

Component	Item Number	Lot Number
SBG Concentrate	100439	821806

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


Review/Approval

Sivasankari Murugan

QC Supervisor

Name

Title

 19 NOV 2018

Signature/Date

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