

Product Number:	102295
Lot Number:	501906
Expiration:	22-Apr-2020

Component	Item Number	Lot Number
SBG Concentrate	103397	911210

SBG Concentrate Concentration (nM)	50
------------------------------------	----

Review/Approval

Patrizia Stadler Assoc. Director, QC

P. Stadler 10 JUL 2019

Name

Title

Signature/Date

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