

Product Number:	102295
Lot Number:	501983
Expiration:	24-Jul-2020

Component	Item Number	Lot Number
SBG Concentrate	103397	920512

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

Review/Approval

Patrizia Stadler	Assoc. Director, QC	
<hr/>	<hr/>	<hr/>
Name	Title	Signature/Date

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