

|                        |             |
|------------------------|-------------|
| <b>Product Number:</b> | 102295      |
| <b>Lot Number:</b>     | 501823      |
| <b>Expiration:</b>     | 15-Apr-2020 |

| <b>Component</b> | <b>Item Number</b> | <b>Lot Number</b> |
|------------------|--------------------|-------------------|
| SBG Concentrate  | 103397             | 910511            |

|                                       |    |
|---------------------------------------|----|
| SBG Concentrate<br>Concentration (nM) | 50 |
|---------------------------------------|----|

Review/Approval

Patrizia Stadler

Assoc. Director, QC

*P. Stadler 07 MAY 2019*

Name

Title

Signature/Date

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