

|                 |             |
|-----------------|-------------|
| Product Number: | 103718      |
| Lot Number:     | 503488      |
| Expiration:     | 11-Jul-2022 |



| Component       | Item Number | Lot Number |
|-----------------|-------------|------------|
| SBG Concentrate | 103650      | 210115     |

|                                    |       |
|------------------------------------|-------|
| SBG Concentrate Concentration (nM) | 134.6 |
|------------------------------------|-------|

| Means of Analysis | Result |
|-------------------|--------|
| HD-1/ HDX         | PASS   |

*Data below represents results generated on the Simoa® HD-1 / HD-X Analyzer.*

Review/Approval

Matt Smart  
Name

QC  
Title

MHA 20 May 2022  
Signature/Date