

| | |
|-----------------|-------------|
| Product Number: | 103718 |
| Lot Number: | 502557 |
| Expiration: | 10-Dec-2020 |

| Component | Item Number | Lot Number |
|-----------------|-------------|------------|
| SBG Concentrate | 103650 | 025414 |

| | |
|------------------------------------|-----|
| SBG Concentrate Concentration (nM) | 135 |
|------------------------------------|-----|

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

Review/Approval

Patrizia Stadler Assoc. Director, QC
Name Title

P. Stadler 01 OCT 2020
Signature/Date