

| | |
|-----------------|-------------|
| Product Number: | 103297 |
| Lot Number: | 135307 |
| Expiration: | 19-Dec-2022 |
| Storage: | 2-8°C |

| Characteristic | |
|------------------------------|------|
| SBG Stock Concentration (nM) | 10nM |

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

Review/Approval

Shivani Goel

QC

Shivani Goel 21 JAN 2022

Name

Title

Signature/Date