

| | |
|------------------------|------------|
| Product Number: | 103397 |
| Lot Number: | 106821 |
| Expiration: | 9-Mar-2022 |
| Storage: | 2-8°C |

Laboratory Analysis

| Characteristic | Result |
|-----------------------|---------------|
| HD-1 Analyzer | PASS |

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

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 13Apr2021
Signature/Date