

| | |
|-----------------|------------|
| Product Number: | 101735 |
| Lot Number: | 502868 |
| Expiration: | 9-Mar-2022 |

| Component | Item Number | Lot Number |
|---------------------|-------------|------------|
| SBG Dilution Buffer | 100376 | 108323 |
| SBG Concentrate | 103397 | 106821 |

Laboratory Analysis

| Means of Analysis | Characteristic | Result |
|------------------------------------|-----------------|--------|
| HD-1 Analyzer | Functional Test | PASS |
| SBG Concentrate Concentration (nM) | | 50 |

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 14Apr2021
Signature/Date