

| | |
|------------------|-----------|
| Product Number: | 101735 |
| Lot Number: | 504215 |
| Expiration Date: | 14-Sep-25 |

| Component | Part Number | Lot Number |
|---------------------|-------------|------------|
| SBG Dilution Buffer | 100376 | 407412 |
| SBG Concentrate | 103397 | 408804 |

Laboratory Analysis

| Characteristic | Means of Analysis | PASS / FAIL |
|----------------|-------------------|-------------|
| Concentration | Spectrophotometer | PASS |

Result

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

Review/Approval

ROBERT SOMOGYI

Name

QC MANAGER

Title

[Signature] 16 APR 2024

Signature/Date