

| | |
|------------------------|-------------|
| Product Number: | 101735 |
| Lot Number: | 502261 |
| Expiration: | 15-Jan-2021 |

| Component | Item Number | Lot Number |
|----------------------------|--------------------|-------------------|
| SBG Dilution Buffer | 100376 | 001516 |
| SBG Concentrate | 103397 | 002714 |

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

Review/Approval Patrizia Stadler Assoc. Director, QC *P. Stadler* 04 Feb 2020
Name Title Signature/Date