

| | |
|------------------------|------------|
| Product Number: | 103397 |
| Lot Number: | 927507 |
| Expiration: | 2-Oct-2020 |
| Storage: | 2-8°C |

Laboratory Analysis

| Characteristic | Result |
|------------------------------------|---------------|
| HD-1 Analyzer | PASS |
| SBG Concentrate Concentration (nM) | 136 |

Review/Approval

Patrizia Stadler

Assoc. Director, QC

P. Stadler

18 Oct 2019

Name

Title

Signature/Date