

Product Number:	103397
Lot Number:	017607
Expiration:	24-Jun-2021
Storage:	2-8°C

Laboratory Analysis

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

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 20 Aug 2020

Name

Title

Signature/Date