

Product Number:	103397
Lot Number:	106103
Expiration:	2-Mar-2022
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
HD-1 Analyzer	PASS

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

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 26 Mar 2021
Signature/Date