

Product Number:	103650
Lot Number:	110217
Expiration:	12-Jul-2021
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
HD-1 Analyzer	PASS

SBG Concentrate Concentration (nM)	152
---------------------------------------	-----

Review/Approval

Shivani Goel

Name

QC

Title

Shivani Goel 07 May 2021

Signature/Date