

Product Number:	103397
Lot Number:	222816
Expiration:	26-Jul-2023
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
HD-1 Analyzer	PASS

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

Review/Approval

Sonia Khillan
Name

QC Manager
Title

Sonia Khillan / 26 OCT 2022
Signature/Date