

Product Number:	103397
Lot Number:	215710
Expiration:	6-Jun-2023
Storage:	2-8°C

Laboratory Analysis

Characteristic	Result
HD-1 Analyzer	PASS

SBG Concentrate Concentration (nM)	58
------------------------------------	----

Review/Approval

Sonia Khilkin
Name

QC Manager
Title

Sonia Khilkin 21MAR2023
Signature/Date