

Product Number:	103397
Lot Number:	530003
Expiration:	27-Apr-2027
Storage:	2-8°C

Laboratory Analysis

Characteristic	Means of Analysis	PASS / FAIL
Functionality	HD(X/1) Analyzer	PASS

RESULT

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

Review/Approval

Kathleen Cornell

Name

Quality Control
Supervisor

Title

[Signature]
20 Nov 2025

Signature/Date