

Product Number:	103397
Lot Number:	534401
Expiration:	10-Jun-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

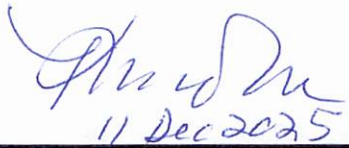
Review/Approval

Kathleen Cornell

Name

Quality Control Supervisor

Title


11 Dec 2025

Signature/Date