

Product Number:	101735
Lot Number:	503788
Expiration:	13-Sep-2023



Data below represents results generated on the Simoa® HD-1 Analyzer.

Component	Item Number	Lot Number
SBG Dilution Buffer	100376	300609
SBG Concentrate	103397	225614

Laboratory Analysis

Means of Analysis	Characteristic	Result
HD-1 Analyzer	Functional Test	PASS
SBG Concentrate Concentration (nM)		50

Review/Approval	<i>Sonia Khillan</i>	<i>QC Manager</i>	<i>Sonia Khillan / 19 JAN 2023</i>
	Name	Title	Signature/Date