

Product Number:	103615
Lot Number:	503758
Paired Reagent Kit	503754
Expiration:	6-Dec-2023
Platform(s):	HD-1 / HD-X & SR-X



Component	Part Number	Lot Number
N2PB Sample Diluent	103516	234010
Calibrator Diluent	103614	234309
NF-light Calibrator Concentrate	102619	220014
GFAP Calibrator Concentrate	103622	219910

Data below represents results generated on the Simoa® HD-1 / HD-X & SR-X Analyzer.

Calibrator Concentrate Value Assignment (pg/mL)	
NF-Light	5240
GFAP	104941

Review/Approval

*Tony Lim*  
Name

QC

Title

*Ronald* 22DEC2022  
Signature/Date