

Product Number:	103671
Lot Number:	503875
Paired Reagent Kit	503819
Expiration:	31-Jan-2024
Platform(s):	HD-1 / HD-X & SR-X



Component	Part Number	Lot Number
N4PE Plasma Sample Diluent	103659	306207
N4PE CSF Sample Diluent	103727	306208
Calibrator Diluent	103661	303111
Abeta 40 Calibrator Concentrate	101734	310203 ¹
Abeta 42 Calibrator Concentrate	103217	211005
GFAP Calibrator Concentrate	103622	219910
NF-light Calibrator Concentrate	102619	220014

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

¹ Control Lot Number for Abeta 40 Calibraor Concentrate - 310203A

Calibrator Concentrate Value Assignment (pg/mL)	
Abeta 40	2478
Abeta 42	2125
GFAP	82774
NF-Light	3580

Review/Approval

Sonia Khillan

Name

QC

Title

Sonia Khillan 11MAY2023

Signature/Date