

Product Number:	103671
Lot Number:	503207
Paired Reagent Kit	503085
Expiration:	26-May-2022
Platform(s):	HD-1 / HD-X & SR-X

Component	Part Number	Lot Number
N4PE Plasma Sample Diluent	103659	121409
N4PE CSF Sample Diluent	103727	114612
Calibrator Diluent	103661	121309
Abeta 40 Calibrator Concentrate	101734	107906
Abeta 42 Calibrator Concentrate	103217	026801
GFAP Calibrator Concentrate	103622	027001
NF-light Calibrator Concentrate	102619	026716

Calibrator Concentrate Value Assignment (pg/mL)	
Abeta 40	4229
Abeta 42	2184
GFAP	149367
NF-Light	9343

Review/Approval

Junpil Park

Name

QC Supervisor

Title

 17 Nov 2021
Signature/Date