

Product Number:	103615
Lot Number:	503172
Paired Reagent Kit	503094
Expiration:	6-Apr-2022
Platform(s):	HD-1 / HD-X & SR-X

Component	Part Number	Lot Number
N2PB Sample Diluent	103516	123011
Calibrator Diluent	103614	109615
NF-light Calibrator Concentrate	102619	026716
GFAP Calibrator Concentrate	103622	027001

Calibrator Concentrate Value Assignment (pg/mL)	
NF-Light	6683
GFAP	133346

Review/Approval

Shivani Goel  
Name

QC Manager  
Title

*Shivani Goel* 04 May 2022  
Signature/Date