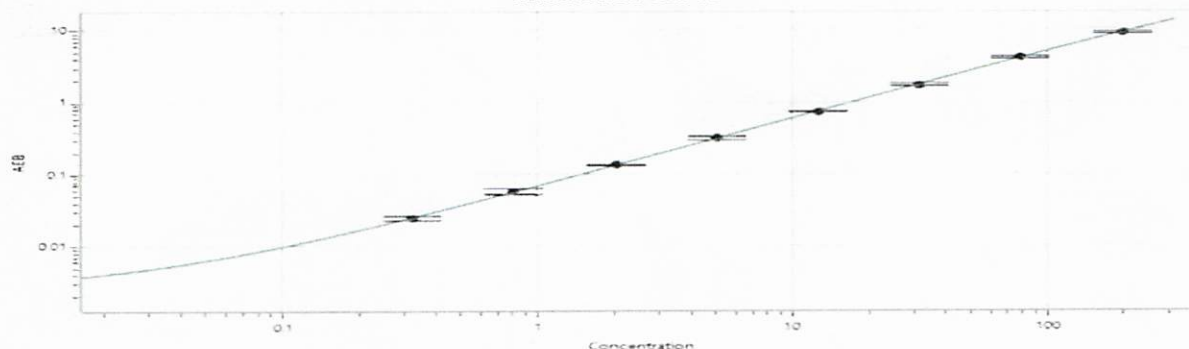


Component	Part Number	Lot Number
Bead Reagent	104778	527504
Detector Reagent	104780	527606
SBG Reagent	104782	527607
Plasma Sample Diluent	104783	527511
CSF Sample Diluent	104784	527515
Calibrator Diluent	104786	527513
Control Diluent	104787	527514
Calibrator Concentrate	104793	409518
Control 1 Concentrate	104795	409519
Control 2 Concentrate	104796	409520
RGP 2.0 Reagent	104321	N/A ¹
Activation Buffer Reagent	104324	N/A ¹

Release Materials*				* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	Spec. Range (pg/mL)	
Control 1 Concentrate	409519	10.8	7.07 - 13.1	
Control 2 Concentrate	409520	459	304 - 564	

**Calibrator Levels
(pg/mL)**

Review/Approval

Dhwani Patel QA Specialist II DP / 19 Nov 2025

Name Title Signature/Date