

Product Number:	104827
Lot Number:	504564
Expiration:	10-Apr-2026

Component	Part Number	Lot Number
Bead Diluent Reagent	104834	410110
Detector Diluent Reagent	104850	412719
SBG Diluent Reagent	104829	410212
Control 1 Diluent Reagent	104841	413409
Control 2 Diluent Reagent	104846	413410
Beads	104833	413809
Detector	104849	413810
SBG	104831	413811
Control 1	104839	408718
Control 2	104844	408719
RGP 2.0 Reagent	104321	N/A ¹
Activation Buffer Reagent	104324	N/A ¹

¹ Reagents are not Kit Lot Specific

Review/Approval

Kyle Giroux

Senior Document Control Specialist, QA

24 Mar 2025

Name

Title

Signature/Date