

Reliable genotyping of APOE-ε4 and other markers in GC-rich regions with a new Axiom Workflow

Monica Chadha, Anu Mittal, Joseph McGirr, Neeraja Vegesna, Carsten Bruckner; Thermo Fisher Scientific, Santa Clara, CA 95051, USA

Abstract

Purpose: Genotyping variants in GC-rich genomic regions remain a challenge for microarrays and short-read sequencing technologies. APOE-ε4 (rs429358), a key genetic risk factor variant for late-onset Alzheimer's disease, has been particularly difficult to genotype due to its location within GC-rich sequences. We have developed a modified whole-genome amplification Axiom assay that not only improves the genotyping performance of APOE but also enhances the detection of other biologically and functionally important variants, especially those located in GC-rich promoters, CpG islands, pharmacogenes, and the MHC- these regions remain technically difficult to genotype with high confidence.

Methods: The study involved amplifying and precipitating gDNA from various sample types using the Axiom™ SwiftArray™ Assay, followed by hybridization and measurement of Axiom™ PangenomePro targets using the SwiftArrayStudio™ Microarray Analyzer. As a representative benchmark, over one hundred SNPs with local GC content ≥ 70% from the 1000 Genomes high coverage call set were evaluated.

Results: The modified workflow achieved a call rate above 99.5%, an average cluster resolution improvement of 1.8 units, and 100% concordance with the 1000 Genomes truth set. Notably, GC-rich loci—including APOE-ε4 (rs429358) and other functionally important variants across key genes such as CYP2A6 (rs1801272, linked to altered nicotine metabolism), KCNQ1 (rs2237897, implicated in type 2 diabetes), and HLA-DPB1 (rs1042151, associated with immune response and aspirin-induced asthma)—showed tighter genotype clusters, improving by 1–4 units compared to the original Axiom protocol. Overall, the modified assay delivered a genome-wide average cluster resolution of ~7.8 units, providing clear evidence that it consistently resolves variants in high-impact genes.

Introduction

The new SwiftArrayStudio Microarray Solution includes the Applied Biosystems™ Axiom™ SwiftArray™ Assay, the Applied Biosystems™ Axiom™ SwiftArrayStudio™ Microarray Analyzer, and the Automated Axiom Analysis Suite as shown in Figure 1.

This advanced solution offers a two-day workflow from sample to insight, featuring improved whole genome amplification that ensures reliable performance for genotyping variants in GC-rich genomic regions while delivering results under 30 hours.

The new assay couples (i) a modified isothermal whole-genome amplification chemistry with enhanced strand-displacement kinetics, (ii) modified enzyme recipe that protects enzyme inhibition and depolymerization via phosphorolysis and (iii) streamlined post-amplification cleanup. The workflow eliminates several Axiom workflow and wash steps, reducing hands-on time to under 3 hours while preserving array throughput. For enhanced sustainability, the solution uses fewer pipette tips with a universal tray pour-and-go mechanism, and the hands-on time is reduced to just 3 hours.

The new workflow exhibits high concordance with the Axiom 2.0 workflow for genotype, copy number call rate, copy number gain/loss, and PGx + other module concordance. Crucially, the new chemistry maintains good representation across the full 20–85 % GC spectrum, yielding tighter genotype clusters.

Materials and Methods

Microarray data: The 96 samples (Blood, Buccal, Saliva, Cell Line) were run on Axiom™ PangenomePro array using Axiom™ SwiftArray™ Assay, Axiom™ SwiftArrayStudio™ Microarray Analyzer and workflow (Figure 1). Sample preparation was done with Axiom™ SwiftArray™ Assay workflow and the current Axiom 2.0 Workflow. All in process QCs met the acceptance criteria. The DAT and CEL files were generated, and Axiom Best Practices Workflow was run to generate genotyping calls. Array genotyping calls were directly compared to 1000 Genomes Phase 3 reference calls to calculate concordance. For matched Blood, Buccal and Saliva samples, genotype call reproducibility was assessed at high GC-rich loci.

Performance evaluation: The performance of genotyping calls from microarray data was assessed for specific markers in GC-rich regions by analyzing the cluster patterns in a two-dimensional space, based on signal intensity and contrast from two channels. Probesets that perform well for a given marker typically exhibit distinct and well-separated clusters. (Figure 2) For rare markers where only a single homozygote cluster is observed, additional samples were examined to confirm that the probeset can make accurate heterozygote calls.

Figure 1. The new Applied Biosystems™ SwiftArrayStudio™ Solution includes:
(1a) Applied Biosystems™ Axiom™ SwiftArray™ Target Preparation method
(1b) Applied Biosystems™ Axiom™ SwiftArrayStudio™ Microarray Analyzer for automated array processing
(1c) Illustrates the advancements of the Axiom™ SwiftArray Assay compared with the Axiom 2.0 Assay.

