

Microarrays

Rapid companion animal genotyping with the SwiftArrayStudio Microarray Analyzer

In this application note, we show:

- A workflow for obtaining companion animal DNA from buccal swabs
- A simplified workflow for collecting genotypes at 630,000 loci of dog and cat genomes
- High-quality genotype information can be obtained from companion animals in less than 30 hours

Introduction

There are an estimated 600 million cats and more than 700 million dogs that are kept as pets worldwide. They are as much a part of our lives as our spouses, our children, and our extended family. We therefore want to understand as much about their health as possible. Advances in genotyping help identify deleterious gene variants that can be important for understanding breeds and for veterinary care.

Genotyping cats and dogs offers significant benefits for both individual animal health and broader population management. For individual owners, genotyping can clarify breed ancestry, informing behavior expectations, nutritional needs, and exercise recommendations [1]. Genetic testing can identify inherited disease risks (e.g., cardiomyopathies, eye diseases, bleeding disorders [1-2]), enabling earlier screening, targeted monitoring, and researching preventive care plans. Genotyping can guide dose selection or drug choice where pharmacogenomic variants are known to affect response to a treatment [3]. For breeders and shelters, genotyping helps reduce the propagation of harmful alleles by informing mating decisions, supporting healthier breeding programs, and giving adopters transparent information about an animal's likely size, temperament, or health trajectory [4-5]. On a population scale, aggregating genomic and phenotype data across millions of pets supports more accurate prevalence estimates for genetic conditions, improves the power of genome-wide association studies, and accelerates the development of new diagnostics and therapeutics in veterinary research [4]. Finally, for public health author-

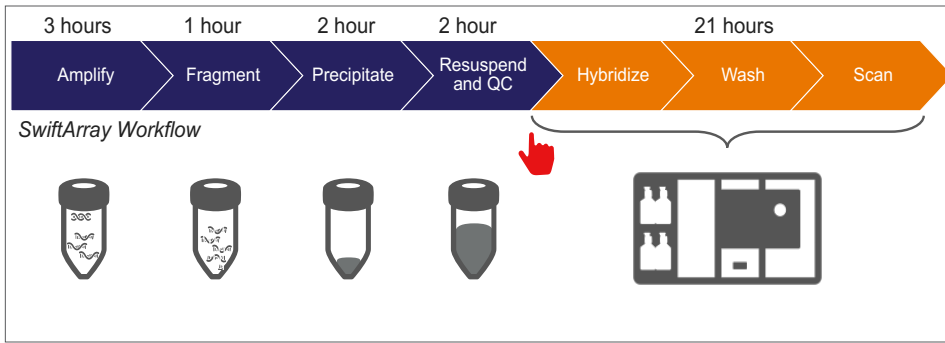


Figure 1. Introducing the SwiftArrayStudio Microarray Analysis solution. The SwiftArray target preparation steps can be completed in about eight hours in a single working day. On-instrument, autonomous array processing can generate data within 21 hours of loading. Overall, a full set of genotypes and copy number calls from an Axiom plate can be generated in about 30 hours. Different colors represent workflow steps on different days; the finger represents a touch point on the instrument, where plates and/or reagents are loaded onto the instrument.

ities, better genetic and demographic data on the companion animal population support more effective vaccination, parasite-control, zoonotic disease surveillance strategies [6-7], and therapeutics in veterinary research [4]. Finally, for public health authorities, better genetic and demographic data on the companion animal population support more effective vaccination, parasite-control, and zoonotic disease surveillance strategies [6-7].

Thermo Fisher Scientific continues to be an innovator of genetic analysis tools. To meet the growing demands for speed in microarray analyses, we recently developed the Applied Biosystems™ SwiftArrayStudio™ Microarray Analyzer and SwiftArray™ Assay. This solution consists of a DNA target preparation workflow that can be completed in a single workday, and a new instrument that automates hybridization, washing, and scanning without any in-process touch points. The SwiftArrayStudio Microarray Analyzer can provide data on an Applied Biosystems™ Axiom™ array plate after an overnight hybridization and scan (Figure 1). This assay and improved workflow reduce the sample-to-answer time from 5 days to 2 days, allowing researchers to gather and interpret genetic data to keep pace with their needs.

Design of experiment

In this application note, we show how the SwiftArrayStudio solution can be used for companion animal genotyping. Briefly, to reflect real-world sample collection conditions, we asked colleagues to obtain buccal samples from their pet cats and dogs using Thermo Scientific™ GenoTube™ Livestock Swabs. Genomic DNA was purified from the swabs using a modified protocol for the Invitrogen™ RecoverAll™ Total Nucleic Acid Isolation kit for FFPE. Up to 100 ng of DNA was processed using the manual SwiftArray Assay workflow in triplicate. Genotypes were obtained using Applied Biosystems™ Axiom™ Feline HD and Axiom™ Canine HD microarray plates. These arrays were designed with extensive consultation with subject matter experts and contain over 630,000 SNP markers that facilitate identification of variants, including those associated with traits like coat color, morphology, and disease risk. For more information, see see our [companion animal web page](#). The prepared samples and plates were loaded onto the SwiftArrayStudio Microarray Analyzer for automated data collection. The data were initially analyzed using Axiom Analysis Suite, and meaningful trait information was obtained from collaborators.

| A. INPUT AMOUNT (ng) | | | | | | | | | | | | |
|----------------------|------|-----|-----|-----|-----|-----|-----|-----|-----|----|-------|----|
| | 1 | 2 | 3 | 4 | 5 | 6 | 7 | 8 | 9 | 10 | 11 | 12 |
| A | >100 | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 70 | 21.4 | 50 |
| B | >100 | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 62 | 19.28 | 25 |
| C | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 44 | 50 | 32 |
| D | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 96 | 44 | 50 | 30 |
| E | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 94 | 44 | 50 | 29 |
| F | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 86 | 43 | 50 | 31 |
| G | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 80 | 42 | 50 | 32 |
| H | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 78 | 32 | 50 | 30 |

| B. Average DNA Amount Recovered (µg) Median = 1067 µg | | | | | | | | | | | | |
|---|--------|--------|--------|--------|--------|--------|--------|--------|--------|--------|--------|--------|
| | 1 | 2 | 3 | 4 | 5 | 6 | 7 | 8 | 9 | 10 | 11 | 12 |
| A | 649.1 | 1101.0 | 1149.3 | 1030.3 | 962.6 | 1144.7 | 1143.3 | 1154.0 | 1119.7 | 1099.7 | 891.1 | 1074.1 |
| B | 812.4 | 1106.7 | 1121.0 | 1089.0 | 1087.7 | 1095.0 | 1108.0 | 1119.0 | 1115.7 | 1069.3 | 825.9 | 1082.7 |
| C | 921.7 | 1064.6 | 1090.1 | 1099.3 | 1034.3 | 1057.9 | 1066.2 | 1060.3 | 1042.9 | 1006.7 | 1060.7 | 962.2 |
| D | 1079.3 | 1039.5 | 1037.7 | 1046.7 | 1012.3 | 1039.7 | 992.3 | 911.5 | 929.9 | 959.8 | 885.0 | 859.9 |
| E | 1127.0 | 1031.5 | 1091.7 | 1072.7 | 1079.3 | 1062.7 | 1051.7 | 1082.0 | 1062.9 | 1041.0 | 770.3 | 1040.0 |
| F | 1080.7 | 1059.6 | 1110.7 | 943.5 | 1087.0 | 1037.6 | 1113.7 | 1083.0 | 1104.7 | 1020.0 | 1117.3 | 1026.8 |
| G | 1140.3 | 1089.2 | 1073.7 | 1093.7 | 1038.9 | 1055.3 | 1105.0 | 1115.7 | 1095.3 | 1099.7 | 1029.8 | 869.1 |
| H | 708.2 | 1016.1 | 1045.8 | 1106.4 | 1079.3 | 1064.1 | 1143.0 | 1126.7 | 1138.3 | 1120.0 | 978.9 | 855.9 |

| C. Average dQC Median = 0.9977 | | | | | | | | | | | | |
|--------------------------------|-------|-------|-------|-------|-------|-------|-------|-------|-------|-------|-------|-------|
| | 1 | 2 | 3 | 4 | 5 | 6 | 7 | 8 | 9 | 10 | 11 | 12 |
| A | 0.551 | 1.000 | 0.982 | 0.999 | 0.582 | 0.995 | 0.999 | 0.998 | 1.000 | 1.000 | 0.998 | 0.998 |
| B | 0.938 | 0.997 | 0.998 | 1.000 | 0.998 | 0.999 | 0.998 | 0.998 | 0.999 | 1.000 | 0.995 | 1.000 |
| C | 0.999 | 0.997 | 0.994 | 0.994 | 0.997 | 0.997 | 0.997 | 1.000 | 1.000 | 1.000 | 1.000 | 0.995 |
| D | 0.988 | 0.993 | 0.996 | 0.999 | 0.998 | 0.992 | 0.999 | 0.999 | 0.999 | 0.996 | 0.997 | 0.787 |
| E | 0.999 | 0.997 | 0.992 | 0.995 | 0.998 | 0.999 | 0.998 | 0.999 | 0.999 | 0.998 | 0.960 | 0.991 |
| F | 0.990 | 0.992 | 0.997 | 0.956 | 0.994 | 0.984 | 0.999 | 0.999 | 1.000 | 0.999 | 0.992 | 0.995 |
| G | 0.998 | 0.985 | 0.993 | 0.995 | 0.995 | 0.993 | 0.999 | 0.989 | 1.000 | 1.000 | 0.999 | 0.322 |
| H | 0.991 | 0.996 | 0.999 | 0.999 | 0.995 | 0.998 | 1.000 | 1.000 | 1.000 | 1.000 | 0.387 | 0.961 |

| D. Average Total Call Rate (%) Median = 99.18 | | | | | | | | | | | | |
|---|--------|--------|--------|--------|--------|--------|--------|--------|--------|--------|--------|--------|
| | 1 | 2 | 3 | 4 | 5 | 6 | 7 | 8 | 9 | 10 | 11 | 12 |
| A | 0.000 | 99.486 | 98.903 | 99.195 | 0.000 | 98.431 | 98.982 | 99.063 | 99.725 | 99.723 | 99.196 | 98.869 |
| B | 0.000 | 99.174 | 99.321 | 99.182 | 99.030 | 99.667 | 97.988 | 99.117 | 99.559 | 99.532 | 98.876 | 99.653 |
| C | 99.233 | 99.221 | 98.531 | 98.423 | 99.004 | 99.460 | 99.024 | 99.368 | 99.603 | 99.270 | 99.515 | 98.347 |
| D | 0.000 | 98.574 | 99.074 | 99.553 | 99.157 | 99.138 | 99.382 | 99.227 | 99.209 | 99.203 | 98.684 | 0.000 |
| E | 99.822 | 99.055 | 98.217 | 99.106 | 99.534 | 99.516 | 98.944 | 98.316 | 98.865 | 99.252 | 0.000 | 0.000 |
| F | 98.346 | 98.333 | 98.757 | 0.000 | 98.225 | 97.505 | 98.950 | 99.512 | 99.443 | 99.354 | 98.903 | 99.433 |
| G | 99.059 | 0.000 | 99.503 | 98.329 | 97.987 | 98.742 | 99.399 | 98.588 | 99.654 | 99.541 | 99.367 | 0.000 |
| H | 99.660 | 98.593 | 99.519 | 98.942 | 99.022 | 98.341 | 99.709 | 99.202 | 99.110 | 99.341 | 0.000 | 0.000 |

Figure 2. Quality control metrics of analyses using Canine Genotyping array. Each sample was analyzed in triplicate on three separate arrays, and the average of the three plates is shown here. (A) Input amount of genomic DNA analyzed (ng). (B) Average amount of DNA recovered from the SwiftArray Assay target preparation workflow. (C) dQC (dish QC, measures signal-to-noise in non polymorphic regions of the genome) of each sample in the three replicates. (D) Average total call rate for each sample in the three replicates.

Recovery of genomic DNA

We collected a variety of representative samples from colleagues and shelters, resulting in 82 dog and 98 cat samples. Buccal swabs were collected from pets and stored at room temperature until used. DNA was recovered manually (i.e., with no automation) using the Ambion™ Recoverall™ FFPE DNA with the following adjustments: 400 µl lysis buffer/sample was used, the swabs were incubated with lysis solution at 65°C for 1 hour, and 200 µl of the resulting lysate was used according to the protocol supplied with the Recoverall kit. Final DNA concentrations were measured with the Invitrogen™ Qubit™ fluorimeter. Recovered amounts ranged from 81 ng to 9.6 µg for cats, and 57.6 ng to 5 µg for dogs.

Target preparation for SwiftArray Assay and loading of SwiftArrayStudio Microarray Analyzer

The SwiftArray assay recommends 100 ng of genomic DNA as an input amount, so samples were diluted to the necessary concentration (5 ng/µl in 20 µl). In cases where not enough DNA was recovered, we used the maximum input volume (20 µl). The DNA quantities input into the SwiftArray Assay are shown in Figures 2 and 3. Note that some of the dog samples (wells C11-H12) were replicates of other samples on the same plate but using less input DNA.

Each sample was analyzed in triplicate on different days using manual SwiftArray workflow as described in the [SwiftArray Assay user guide](#). As described in the user guide, it is recommended that the amount of DNA produced and recovered be quantified before proceeding with the hybridization. The average amount of DNA recovered from the dog and cat SwiftArray assay triplicates is shown. Plates are recommended to have a median recovery of around 1000 µg/well; the dog plates had average three-plate median of 1067 µg and the cat plates had average three-plate median of 1066 µg. Furthermore, analysis of fragment sizes by agarose gel electrophoresis showed the samples had correct size range of DNA fragments (data not shown).

Having established that the samples met quality standards, they were prepared and loaded on the SwiftArrayStudio Microarray Analyzer for automated hybridization (16 hour hybridization time), washing and scanning. The total time on the instrument was about 21 hours.

| A. INPUT AMOUNT (ng) | | | | | | | | | | | | |
|----------------------|-----|-----|-----|-----|-----|-----|-----|-----|-----|-----|------|------|
| | 1 | 2 | 3 | 4 | 5 | 6 | 7 | 8 | 9 | 10 | 11 | 12 |
| A | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 91.6 | 71 |
| B | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 90.8 | 52.8 |
| C | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 86.2 | 47.2 |
| D | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 85.6 | 45.6 |
| E | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 80.4 | 43.8 |
| F | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 88.8 | 39.8 |
| G | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 82.0 | 77.2 |
| H | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 100 | 91.8 | 76.2 |
| | | | | | | | | | | | | 27 |

| B. Average DNA Amount Recovered (µg) Median = 1067 µg | | | | | | | | | | | | |
|---|--------|--------|--------|--------|--------|--------|--------|--------|--------|--------|--------|--------|
| | 1 | 2 | 3 | 4 | 5 | 6 | 7 | 8 | 9 | 10 | 11 | 12 |
| A | 920.5 | 717.0 | 761.7 | 880.2 | 852.3 | 727.2 | 801.8 | 852.4 | 793.5 | 676.8 | 895.9 | 588.1 |
| B | 1062.9 | 1101.5 | 1164.0 | 1080.3 | 1156.7 | 1154.7 | 1165.7 | 1168.3 | 1170.0 | 1173.3 | 1173.0 | 1136.7 |
| C | 854.8 | 817.5 | 890.8 | 963.5 | 961.4 | 956.6 | 956.9 | 948.2 | 961.9 | 906.3 | 898.4 | 934.6 |
| D | 816.9 | 886.2 | 1173.9 | 1055.3 | 1143.0 | 1084.0 | 1092.0 | 1153.3 | 1153.7 | 1182.0 | 1159.7 | 1041.9 |
| E | 955.0 | 1097.3 | 1107.3 | 1125.0 | 1150.3 | 1051.7 | 1116.0 | 1122.7 | 1062.3 | 1069.3 | 1002.0 | 1064.0 |
| F | 816.6 | 1188.3 | 1222.0 | 1161.3 | 1176.3 | 1172.7 | 1173.7 | 1036.1 | 1142.3 | 1202.7 | 1186.7 | 1236.7 |
| G | 794.2 | 1141.3 | 1148.7 | 1154.7 | 1137.3 | 1134.0 | 1086.3 | 1126.7 | 1125.3 | 1146.7 | 1100.3 | 1155.0 |
| H | 635.4 | 850.1 | 1033.3 | 868.5 | 920.5 | 951.9 | 946.5 | 1243.0 | 942.5 | 997.3 | 722.5 | 722.1 |

| C. Average DQC Median = 0.9963 | | | | | | | | | | | | |
|--------------------------------|--------|--------|--------|--------|--------|--------|--------|--------|--------|--------|--------|--------|
| | 1 | 2 | 3 | 4 | 5 | 6 | 7 | 8 | 9 | 10 | 11 | 12 |
| A | 0.9630 | 0.9667 | 0.9927 | 0.9810 | 0.9913 | 0.9963 | 0.9917 | 0.9227 | 0.9823 | 0.9790 | 0.9980 | 0.8370 |
| B | 0.9980 | 0.9897 | 0.9977 | 0.9893 | 0.9973 | 0.9943 | 0.9927 | 0.9973 | 0.9973 | 0.9990 | 0.9980 | 0.9990 |
| C | 0.8017 | 0.9960 | 0.9703 | 0.9937 | 0.9977 | 0.9937 | 0.9960 | 0.9983 | 0.9957 | 0.9980 | 0.9997 | 0.9983 |
| D | 0.9890 | 0.9857 | 0.9967 | 0.9970 | 0.9963 | 0.9947 | 0.9987 | 0.9977 | 1.0000 | 0.9980 | 0.9980 | 0.9990 |
| E | 0.0360 | 0.9967 | 0.9733 | 0.9837 | 0.9973 | 0.9913 | 0.9970 | 0.9963 | 0.9990 | 1.0000 | 0.9990 | 0.9980 |
| F | 0.9825 | 0.9903 | 0.9967 | 0.9970 | 0.9980 | 0.9957 | 1.0000 | 0.5820 | 0.9980 | 0.9993 | 1.0000 | 0.9983 |
| G | 0.9850 | 0.9947 | 0.9927 | 0.9990 | 0.9973 | 0.9987 | 1.0000 | 0.9973 | 0.9990 | 0.9993 | 0.9913 | 0.9960 |
| H | 0.9223 | 0.9957 | 0.9930 | 0.9950 | 0.9970 | 0.9967 | 0.9910 | 0.9960 | 0.9903 | 0.9970 | 0.9950 | 0.0360 |

| D. Average Total Call Rate (%) Median = 99.70 | | | | | | | | | | | | |
|---|--------|--------|--------|--------|--------|--------|--------|--------|--------|--------|--------|--------|
| | 1 | 2 | 3 | 4 | 5 | 6 | 7 | 8 | 9 | 10 | 11 | 12 |
| A | 99.882 | 98.718 | 98.530 | 0.000 | 99.247 | 99.671 | 99.310 | 97.687 | 98.454 | 97.987 | 99.736 | 98.638 |
| B | 99.916 | 99.113 | 99.785 | 98.620 | 99.435 | 99.597 | 99.282 | 99.874 | 99.806 | 99.887 | 99.836 | 99.852 |
| C | 99.902 | 99.726 | 98.695 | 99.385 | 99.837 | 99.465 | 99.767 | 99.789 | 99.347 | 99.821 | 99.807 | 99.874 |
| D | 99.922 | 98.845 | 99.881 | 99.857 | 99.487 | 99.397 | 99.878 | 99.543 | 99.892 | 99.749 | 99.915 | 99.801 |
| E | 0.000 | 99.781 | 97.659 | 97.327 | 99.823 | 99.080 | 99.566 | 99.446 | 99.894 | 99.907 | 99.512 | 99.896 |
| F | 97.545 | 99.082 | 99.748 | 99.908 | 99.809 | 99.660 | 99.884 | 0.000 | 99.798 | 99.913 | 99.907 | 99.860 |
| G | 98.535 | 99.663 | 99.342 | 99.887 | 99.913 | 99.866 | 99.903 | 99.869 | 99.878 | 99.816 | 99.305 | 99.651 |
| H | 98.854 | 99.634 | 99.370 | 99.581 | 99.802 | 99.648 | 99.404 | 99.673 | 99.268 | 99.559 | 99.487 | 0.000 |

Figure 3. Quality control metrics of analyses using the Feline Genotyping array. Each sample was analyzed in triplicate on three separate arrays, and the average of the three plates is shown here. **(A)** Input amount of genomic DNA analyzed (ng). **(B)** Average amount of DNA recovered from the SwiftArray Assay target preparation workflow. **(C)** Average dQC (dish QC, measures signal-to-noise in non polymorphic regions of the genome) of each sample in the three replicates. **(D)** Average total call rate for each sample in the three replicates.

Scanning quality metrics

Following hybridization, washing and scanning on the SwiftArrayStudio Microarray Analyzer, we analyzed the QC metrics generated by the Applied Biosystems™ Axiom™ Analysis Suite and default best practices parameters (see the [Axiom Analysis Suite User Guide](#)). One of these, DishQC (DQC), measures the amount of overlap between two homozygous peaks created by non polymorphic probes, with a value of 1.0 considered optimal. Among the canine samples, six failed the default DQC metric (0.82), corresponding to an approximate 93.8% pass rate. For the cat samples, four failed to meet the default DQC metric (passing rate 95.8%).

Another quality measurement generated by Axiom Analysis Suite is the average passing QC call rate (average percentage of autosomal SNPs with a call other than NoCall, as measured at the sample QC step of passing samples within a plate). Among the dog samples that passed DQC, six more failed the default call rate threshold (97%); similarly, one additional cat sample failed to meet the threshold. However, in the other samples the data quality was high, giving median total call rates of 99.18% for passing dog samples and 99.67% for passing cat samples. Notably, these passing samples included several with substantially less than the recommended 100 ng DNA input.

Table 1. Performance of SwiftArray solution for canine and feline microarrays. Values shown are the average of three triplicate plates analyzing buccal swabs of pet DNA. Abbreviations: DQC: Dish QC; CR: call rate. For more details, see the [Axiom Analysis Suite User Guide](#).

| Metric | Cat Samples | Dog Samples |
|---|-------------|-------------|
| Total number of markers | 662,069 | 712,332 |
| Samples with DQC \geq 0.82 | 95.8% | 93.8% |
| DQC median | 0.9963 | 0.9977 |
| Median passing QC CR | 99.7% | 98.2% |
| Average number of samples with passing QC CR | 95.8% | 87.5% |
| Percent of markers with high performance in dataset | 97.2% | 93.8% |
| Average sample reproducibility | 99.99% | 99.93% |
| Gender calls (all three replicates) | | |
| Female | 136 | 110 |
| Male | 118 | 120 |
| Axiom Analysis Suite displayed values | | |
| PolyHighResolution | 89.29% | 69.83% |
| NoMinorHom | 5.59% | 12.79% |
| MonoHighResolution | 2.30% | 11.15% |

Table 2. Top 20 dog breeds represented in a collection of pets.

Note that most of these are due to mixtures of characteristic loci detected that were in the dogs, not necessarily pure breeds.

| Breed | Number |
|--------------------------------|--------|
| Toy Poodle | 12 |
| Golden Retriever | 11 |
| Vizsla | 9 |
| American Cocker Spaniel | 8 |
| East European Shepherd | 8 |
| American Staffordshire Terrier | 7 |
| Chihuahua | 7 |
| German Shepherd | 7 |
| Laborador Retriever | 7 |
| Poodle | 7 |
| English Shepherd | 5 |
| Siberian Husky | 5 |
| Staffordshire Bull Terrier | 5 |
| Akbash | 4 |
| Central Asian Shepherd Dog | 4 |
| Italian Greyhound | 4 |
| Australian Cattle Dog | 3 |
| Blue Gascony Griffon | 3 |
| Bull Terrier | 3 |
| English Cocker Spaniel | 3 |

Assessment of genotypic clustering quality showed that 91.5% of canine markers were classified as polymorphic high resolution, no minor homozygotes, or monoallelic high resolution (Table 1), while 95.4% of feline markers fell into these clustering categories. For both sets of data, the remainder fell into less informative classes of genotyping clusters (CallRateBelowThreshold, OffTargetVariant, OtherMA or Other). Thus, despite the variable quality of these samples, the SwiftArray assay workflow can generate high-quality genotyping data.

Converting genotypes to traits

The goal of obtaining genotyping data is to gain information on traits, breeds, pathologies, etc. To convert the genotype data to possible phenotypes, we relied on collaborators. For dogs, we sent the data to a collaborator who had developed a custom breed identification pipeline, matching genotypes at specific loci to over 200 dog breeds. Using the data collected from colleagues' dogs, we tallied the most common breed loci identified in these samples (Table 2). Note these do not represent the number of pure-breed dogs; rather, the list indicates which breeds are likely to have contributed to the makeup of the mixed-breed pets owned by our colleagues.

Similarly, an external collaborator provided a more detailed conversion of selected genotypes to selected cat traits (Table 3). Included in this data set were alleles for known genetic pathologies (not shown); fortunately, none of the cats were positive for these alleles. Note that for both the canine and feline arrays, more than 630,000 markers were analyzed; the results shown here are a small fraction of the total collected. For more information on the loci contained on the feline and canine microarrays, see our [companion animal web page](#).

Table 3. Tally of selected cat traits in a collection of 92 cats. Some traits and genes are listed twice; this due to the two different alleles of the same gene being examined. Note that this is a subset of the more than 640,000 genotypes analyzed on the array.

| Trait | Gene | # Ref homo | # Ref hetero | # Alt homo |
|--------------------------------|--------------|------------|--------------|------------|
| Burmese coat color | <i>TRY</i> | 90 | 1 | 1 |
| Ticked | <i>DKK4</i> | 85 | 4 | 3 |
| Polydactyly | <i>LMBR1</i> | 90 | 2 | 0 |
| Sphynx Hairless | <i>KRT71</i> | 70 | 19 | 1 |
| Chocolate | <i>TYRP1</i> | 85 | 7 | 0 |
| Coat color, dilute | <i>MLPH</i> | 35 | 33 | 24 |
| Long hair | <i>FGF5</i> | 64 | 25 | 3 |
| Long hair | <i>FGF5</i> | 49 | 32 | 11 |
| Blotched tabby | <i>LVRN</i> | 22 | 41 | 28 |
| Coat color, non-agouti (black) | <i>ASIP</i> | 8 | 35 | 49 |
| Coat color, Salmiak | <i>KIT</i> | 92 | 0 | 0 |
| Curly coat, Devon Rex | <i>KRT71</i> | 66 | 12 | 10 |
| Curly coat, Ural Rex | <i>LIPH</i> | 92 | 0 | 0 |
| Coat color, russet | <i>MC1R</i> | 92 | 0 | 0 |
| Coat color, copal | <i>MC1R</i> | 92 | 0 | 0 |
| Tail, short | <i>TBXT</i> | 92 | 0 | 0 |
| Tail, short | <i>TBXT</i> | 91 | 1 | 1 |
| Munchkin, standard | <i>UGDH</i> | 92 | 0 | 0 |

Summary and conclusions

In this application note, we describe a simplified method for obtaining genotypic information from dogs and cats. We showed how genomic DNA can be collected from buccal swabs using GenoTube swabs, and purified using RecoverAll kits. We also showed how the SwiftArrayStudio Microarray Analyzer and Assay can be used to generate high quality genotypic data from a variety of real-world users. Importantly, the SwiftArray Assay can generate these data in less than 30 hours. The SwiftArray-Studio Microarray Processor Solution, together with Axiom Feline and Canine microarrays, is therefore poised to provide breeders, veterinarian researchers, and providers of genetic data with tools that can generate high-quality genotyping information rapidly, accurately, and reliably.

To obtain the library files optimized for the SwiftArrayStudio Microarray Analyzer, please speak with your local FAS.

Acknowledgements

We thank Dr. Leslie Lyons at the University of Missouri for her assistance in designing the feline array and for analyzing the pet genotypes generated in this study. We thank Dr. Dongyeop Kim for his analysis of the dog genotypes presented here.

References

1. Lyons LA. Genetic Testing: practical dos and don'ts for cats. *J Feline Med Surg.* 2024 Dec;26(12):1098612X241303603. doi: 10.1177/1098612X241303603.
2. Dreger DL, Rimbault M, Davis BW, Bhatnagar A, Parker HG, Ostrander EA. Whole-genome sequence, SNP chips and pedigree structure: building demographic profiles in domestic dog breeds to optimize genetic-trait mapping. *Dis Model Mech.* 2016 Dec 1;9(12):1445-1460. doi: 10.1242/dmm.027037.
3. Champion DP, Dowell FJ. Translating Pharmacogenetics and Pharmacogenomics to the Clinic: Progress in Human and Veterinary Medicine. *Front Vet Sci.* 2019 Feb 11;6:22. doi: 10.3389/fvets.2019.00022.
4. Mousa Basha SA, Akis I. Missense variants in MC4R gene are associated with obesity in cats. *Vet Res Commun.* 2025 Mar 4;49(3):127. doi: 10.1007/s11259-025-10700-4.
5. Ludwig-Peisker O, Ansel E, Schweizer D, Jagannathan V, Loechel R, Leeb T. PCYT1A Missense Variant in Vizslas with Disproportionate Dwarfism. *Genes (Basel).* 2022 Dec 13;13(12):2354. doi: 10.3390/genes13122354.

Ordering information

| Description | Cat. No |
|--|---------|
| SwiftArrayStudio™ Microarray Analyzer | 00-0471 |
| Axiom™ SwiftArrayStudio Consumables Kit | 952674 |
| Axiom™ Feline HD Array | 551642 |
| Axiom™ Canine HD Array | 550869 |
| Thermo Scientific™ GenoTube™ Livestock Swabs | 9062010 |
| Invitrogen™ RecoverAll™ Total Nucleic Acid Isolation Kit for FFPE | AM1975 |
| Invitrogen™ E-Gel™ 48 Agarose Gel with SYBR Safe stain, 4% agarose | G820804 |
| 25-bp DNA Ladder | 931343 |

 Learn more at thermofisher.com/companionanimal

For Research Use Only. Not for use in diagnostic procedures. © 2026 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. APN-14327008 0526

applied biosystems