

Attend these presentations during the AACR Annual Meeting 2022 to see how NGS is advancing oncology research

Detection of gene fusions and exon skipping events in lung FFPE samples with the OncoPrint™ Precision Assay on the Ion Torrent™ Genexus™ System

Number: 78 | Board 15

Presenter: Amir Marcovitz

Session title: Cancer Genomics 2

Time and location: Sunday, April 10, 2022
1:30 p.m.–5:00 p.m., Hall D-H, Poster Section 4
Morial Convention Center

High-throughput NGS for the study of oncology variants, gene fusion events, and key oncology research endpoints

Number: 42 | Board 3

Presenter: Dinesh Cyanam

Session title: Cancer Genomics 1

Time and location: Sunday, April 10, 2022
1:30 p.m.–5:00 p.m., Hall D-H, Poster Section 3
Morial Convention Center

Analytical performance of a novel next-generation sequencing assay for myeloid MRD research

Number: 759 | Board 15

Presenter: Wally Zhang

Session title: Cancer Genomics 3

Time and location: Monday, April 11, 2022
9:00 a.m.–12:30 p.m., Hall D-H, Poster Section 3
Morial Convention Center

Fully automated sample-to-report NGS workflow for comprehensive genomic profiling of myeloid neoplasms

Number: 764 | Board 20

Presenter: Sarah Brozio

Session title: Cancer Genomics 3

Time and location: Monday, April 11, 2022
9:00 a.m.–12:30 p.m., Hall D-H, Poster Section 3
Morial Convention Center

Shannon entropy of mutational signatures predicts sensitivity of signature detection in targeted sequencing

Number: 1515 | Board 16

Presenter: Fiona Hyland

Session title: Genomic Instability 2

Time and location: Monday, April 11, 2022
1:30 p.m.–5:00 p.m., Hall D-H, Poster Section 5
Morial Convention Center

Evaluation of multiple myeloma research samples by analysis of multiple B cell receptor targets in a single NGS assay

Number: 2293 | Board 24

Presenter: Geoff Lowman

Session title: Cancer Genomics 4

Time and location: Tuesday, April 12, 2022
9:00 a.m.–12:30 p.m., Hall D-H, Poster Section 5
Morial Convention Center

Automation meets reliability: use of the OncoPrint Precision Assay on the Genexus System for accurate identification of cancer biomarkers in FFPE and liquid biopsy samples

Number: 2939 | Board 11

Presenter: Thilanka Jayaweera

Session title: Cancer Genomics 5

Time and location: Tuesday, April 12, 2022
1:30 p.m.–5:00 p.m., Hall D-H, Poster Section 2
Morial Convention Center

Somatic hypermutation analysis of chronic lymphocytic leukemia research samples by DNA or RNA input IGH chain sequencing

Number: 2940 | Board 12

Presenter: Michelle Toro

Session title: Cancer Genomics 5

Time and location: Tuesday, April 12, 2022
1:30 p.m.–5:00 p.m., Hall D-H, Poster Section 2
Morial Convention Center

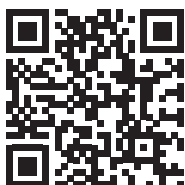
Rapid and accurate variant calling of FFPE samples with the Genexus System

Number: 2944 | Board 16

Presenter: Iris Casuga

Session title: Cancer Genomics 5

Time and location: Tuesday, April 12, 2022
1:30 p.m.–5:00 p.m., Hall D-H, Poster Section 2
Morial Convention Center



Learn more at thermofisher.com/aacr