

## Poster Index:

- Poster 1** Kwasi Agbleke  
**Live Cell Imaging of Homologous Chromosomes During Differentiation**
- Poster 2** Ravindra Amunugama  
**The Role of Replication Fork Regression During Incision-dependent DNA Interstrand Crosslink Repair**
- Poster 3** Maria Andrianova  
**Changes in mutational processes and patterns of mutations during Cancer Development**
- Poster 4** Artem Artemov  
**Cancer genomes revealed the role of POLD in replication of both the leading and the lagging strands at origins**
- Poster 5** Elizabeth Boehm  
**Ubiquitin regulates the catalytic activity of DNA polymerase eta through a novel binding site on the catalytic core**
- Poster 6** Simon Brunner  
**Reconstructing breakpoints of cancer genomes at base-pair resolution.**
- Poster 7** Chongyi Chen  
**Single-Cell Whole Genome Analyses by Linear Amplification via Transposon Insertion (LIANTI)**
- Poster 8** Jia-Yun Chen  
**Molecular dynamics of oncogenic BRaf-induced senescence**
- Poster 9** Lin Deng  
**Complex chromosomal rearrangements from replication errors in mitosis**
- Poster 10** Alan Jiao  
**NRDE2 is an essential gene required for genome stability in human cells**
- Poster 11** Mitchell Leibowitz  
**Defining the genomic consequences and protective mechanisms underlying DNA replication stress**
- Poster 12** Katherine Marcelain  
**Defective H3K9 trimethylation of pericentromeric satellite regions leads to aneuploidy in Ski-deficient cells.**
- Poster 13** Isabella Marcomini  
**Understanding the determinants of telomeric identity for repair pathways**
- Poster 14** So Hae (Irene) Park  
**Role of transcription in non-recurrent copy number variant (CNV) formation**

- Poster 15**                   Dmitriy Podolskiy  
**Analysis of cancer genomes reveals basic features of human aging and its role in cancer development**
- Poster 16**                   Steven Roberts  
**Mechanisms that establish an APOBEC mutation signature in cancer and yeast**
- Poster 17**                   Srinivas Vinod Saladi  
**Transcriptional reprogramming by ACTL6A, SWI/SNF complex subunit, drives YAP activation and regenerative proliferation in HNSCC**
- Poster 18**                   Dan Semlow  
**NEIL3-mediated DNA interstrand cross-link repair is functionally coupled to the replicative helicase**
- Poster 19**                   Maxwell Sherman  
**How do we solve it: PCR Amplification Confounds Copy Number Variation Detection in Whole Genome Sequencing**
- Poster 20**                   Ofar Shoshani  
**Functional Chromothripsis Drives Drug Resistance in Cancer Cells**
- Poster 21**                   Mar Soto Ruiz de la Torre  
**Tolerance to whole chromosome versus partial aneuploidies is determined by their differential effect on genomic stability**
- Poster 22**                   Neil Umbreit  
**Chromosome bridge resolution requires mechanical forces from actin-based contractility**
- Poster 23**                   Jeroen van den Berg  
**A Cas9-based system to study location-dependent effects on the DNA damage response**
- Poster 24**                   Peter Waltman  
**The role and signature of chromosomal instability in Fanconi Anemia Head and Neck Squamous Carcinoma**
- Poster 25**                   Nicholas Willis  
**BRCA1 suppresses tandem duplications at Tus/Ter-stalled replication forks in mammalian cells.**
- Poster 26**                   Hong Xu  
**Targeting DNA G-quadruplex sequences for selective toxicity to BRCA1/2 deficient tumors**
- Poster 27**                   Xiaotong Yao  
**Gnome/SkyWalker suite represents rearranged cancer genome as graph and reconstructs local karyotypes**
- Poster 28**                   Vladimir seplyarskiy  
**Error-prone damage bypass during lagging strand replication is a common source of mutations in cancers and germline**
- Poster 29**                   Ruth Brann McCole  
**Ultraconservation of DNA sequence provides a new lens for focusing on chromosomal structural rearrangements in neurodevelopmental disorder genomes**