



# Whole Genome Sequencing as a Screening Tool in Healthy Population: Lesson learned from 110 cases

Madhuri Hegde, PhD., FACMG, Ruby Liu, Christin Collins, Babi Nallamilli, Megan Landsverk, Jing Xie, Akanchha Kesari, Alice Tanner, Zeq Ma, and Alka Chaubey.  
Revvity Omics, Pittsburgh, PA

## BACKGROUND

- Whole genome sequencing (WGS) has been widely considered a powerful tool in identification of disease-causing variants to ascertain clinical diagnoses.
- With WGS's uniformed coverage for both exonic and intronic regions, the expected clinical sensitivity of WGS is up to 40-50%.
- the utilization of WGS as a screening test for healthy individuals has increased.
- RVTY healthy WGS includes copy number variation analysis and mitochondrial DNA analysis.
- We have performed 110 WGS on reportedly healthy asymptomatic individuals

## RESULTS

Figure 1: Healthy WGS overall yield

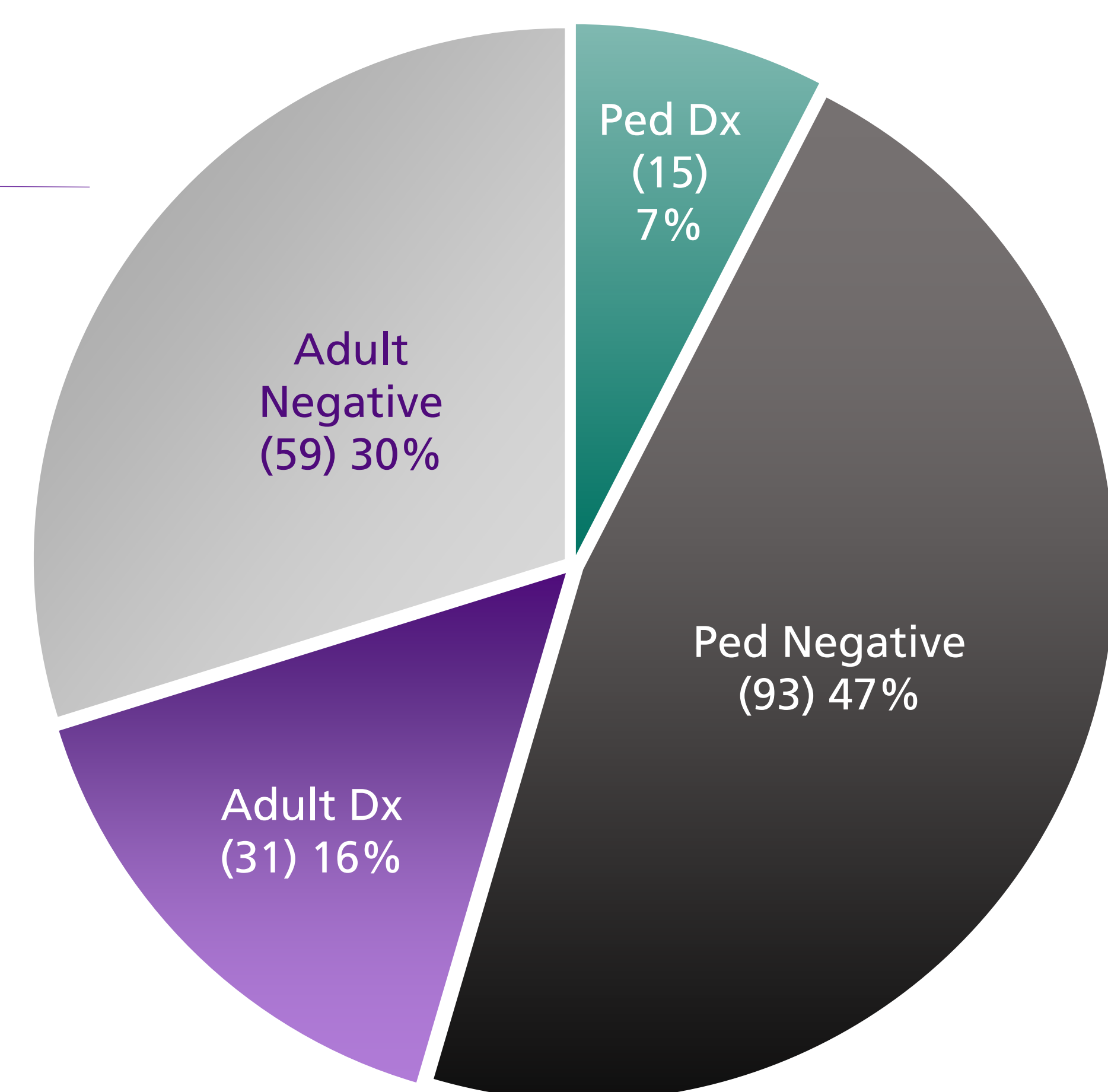


Figure 2: Type of variants in Pediatric and Adult Dx Findings

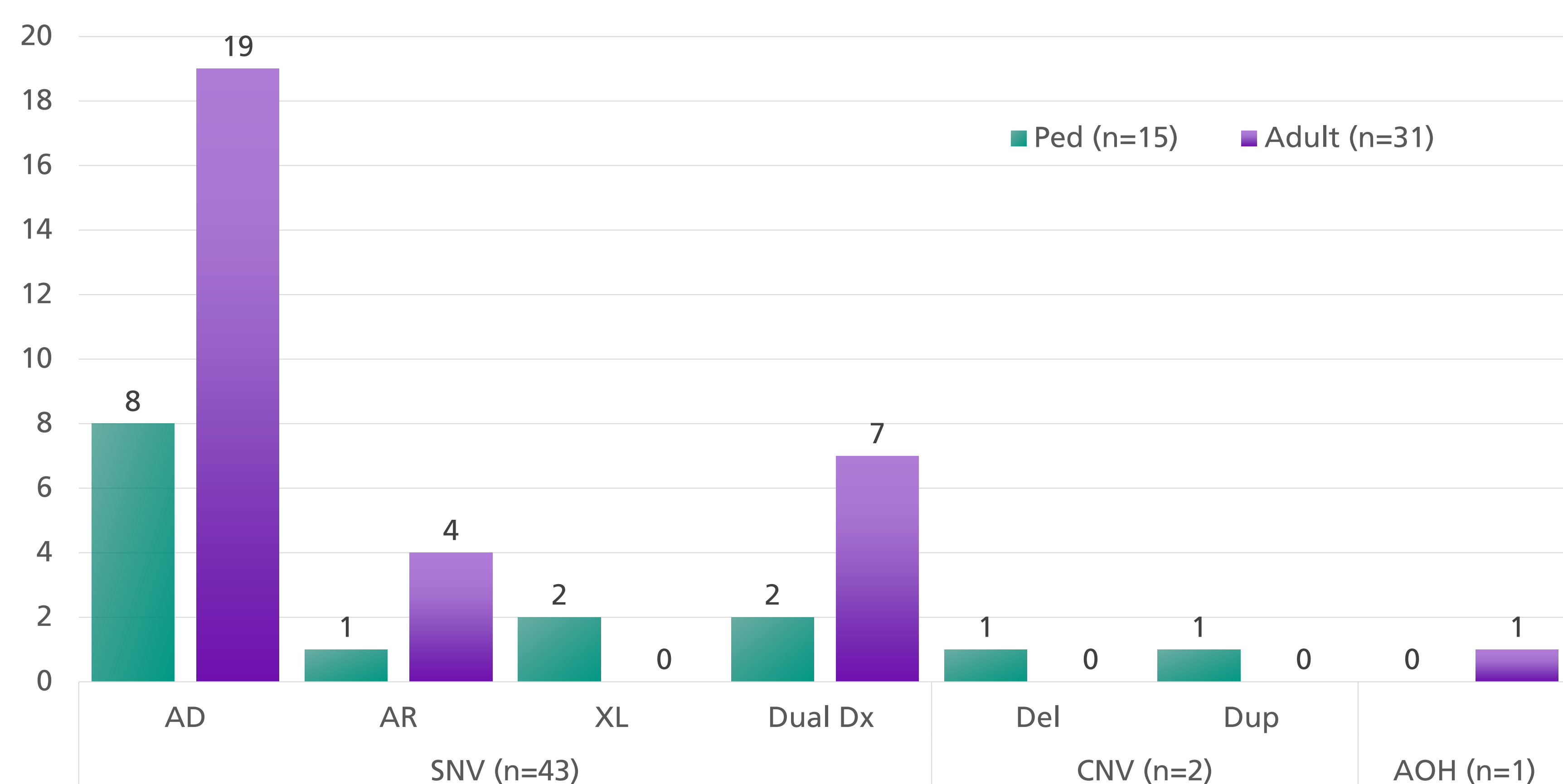


Figure 3: Associated Diseases in Pediatric and Adult Cohort

Category	# of genes	Genes	# of cases
Hereditary cancer syndrome	7	<i>SDHA, MTF, CHEK2, RAD50, FH, APC, PTEN</i>	12
Neurodegeneration	2	<i>APOE, LRRK2</i>	11
Neurology	2	<i>PRRT2, MME</i>	2
Endocrinology	3	<i>ABCC8, PROKR2, NOBOX</i>	3
Skeletal dysplasia	2	<i>ARSE, FGFR3</i>	2
Biochemical	3	<i>HFE, GALT, G6PD</i>	7
Pulmonary	2	<i>SFTPC, SMAD9</i>	2
Urology	2	<i>SLC34A1, SLC3A1</i>	2
Cardiovascular	3	<i>MFAP5, MYBPC3, PRKAG2</i>	4
Muscular	2	<i>AMPD1, SCN4A</i>	3
Other (deafness, skin, dysmorphism, fever)	4	<i>IRF6, FLG, TECTA, MEFV</i>	4

## MULTIPLE DIAGNOSES EXAMPLES

Case/Age	Dx findings 1	Associated disease	Dx findings 2	Associated disease
1/PED	<i>AMPD1</i> , c.[133C>T;242C>T] (p.[Gln45Ter; Pro81Leu]), Pathogenic with reduced penetrance, homozygous	Myopathy due to myoadenylate deaminase deficiency	<i>SDHA</i> , c.1534C>T (p.Arg512Ter), Likely Pathogenic	Parangangliomas 5
2/PED	<i>HFE</i> , c.187C>G (p.His63Asp); c.845G>A (p.Cys282Tyr), Pathogenic with reduced penetrance	Hemochromatosis	<i>TECTA</i> , c.2342_2345dup, Pathogenic	Deafness, autosomal dominant 8/12
3/ADULT	<i>LRRK2</i> , c.6055G>A(p.Gly2019Ser), Pathogenic	Parkinson disease	<i>APOE</i> , c.388T>C (p.Cys130Arg), risk factor	Alzheimer disease
4/ADULT	<i>MYBPC3</i> , c.3697C>T (p.Gln1233Ter), Pathogenic.	Hypertrophic cardiomyopathy; Dilated cardiomyopathy; Left ventricular noncompaction 10	<i>SLC3A1</i> , c.1354C>T (p.Arg452Trp), Pathogenic	Cystinuria
4/ADULT	<i>MME</i> , c.467del, Pathogenic	Charcot-Marie-Tooth disease, axonal, type 2T; Spinocerebellar ataxia 43	<i>PROKR2</i> , c.254G>A (p.Arg85His), Pathogenic	Hypogonadotropic hypogonadism 3 with or without anosmia
5/ADULT	<i>APC</i> , c.3920T>A (p.Ile1307Lys), Risk Factor.	Familial adenomatous polyposis (FAP)	<i>G6PD</i> , c.563C>T (p.Ser188Phe), Pathogenic, Hemizygous	Hemolytic anemia, G6PD deficient (favism)
6/ADULT	<i>MYBPC3</i> , c.3190+5G>MYBPC3 A, Pathogenic	Hypertrophic cardiomyopathy; Dilated cardiomyopathy; Left ventricular noncompaction 10	<i>NOBOX</i> , c.1295_1310delinsGAG, Likely Pathogenic	Premature ovarian failure 5
7/ADULT	<i>PTEN</i> c.801+1del, Pathogenic	Cowden syndrome 1; Macrocephaly/autism syndrome; Lhermitte-Duclos syndrome	<i>PRKAG2</i> c.905G>A(p.Arg302Gln), Pathogenic	Cardiomyopathy, hypertrophic 6; Glycogen storage disease of heart, lethal congenital; Wolff-Parkinson-White syndrome
8/ADULT	<i>MITF</i> , c.952G>A (p.Glu318Lys), Pathogenic	Melanoma, cutaneous malignant, susceptibility to, 8	<i>APOE</i> , c.388T>C (p.Cys130Arg), risk factor	Alzheimer disease

## CONCLUSION

- Healthy WGS screening test can be considered as a new standard of care along with newborn screening and carrier screening.
- Healthy WGS allows individuals to learn their risk for medically actionable conditions and carrier status prior to family planning
- Healthy WGS provide valuable information for the healthcare providers to help their patients making suitable prophylactic plans, finding medical management guidelines, selecting optimal dosage of medications and adapting to a healthy life style.