

Diagnostic Yield and Clinical Utility of Nephrolithiasis and Primary Hyperoxaluria Sequencing panels

Christin D Collins, PhD, FACMG; Yang Wang, PhD, FACMG; Michelle D. Glover, PhD; Madhuri Hegde, PhD, FACMG

revvity



Dicerna

Introduction

- Nephrolithiasis (so called kidney stones, renal calculi or urolithiasis) is one of the most common renal complications with more than 200,000 cases diagnosed in the US every year.
- The etiology of nephrolithiasis is multifactorial involving both genetic and environmental factors.
- Half of all affected individuals have a family history of kidney stones suggesting the genetic trait plays a vital role in disease manifestation.
- PKIG designed a 35-gene Nephrolithiasis and a 3-gene Primary Hyperoxaluria gene panels to identify the genetic underlining of these diseases

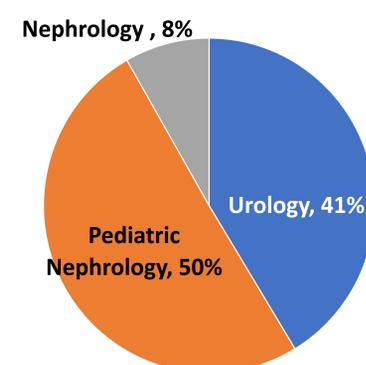
RESULTS

- A diagnosis was reached in 15% of these cases (24 cases) with equal split between Nephrolithiasis and Primary Hyperoxaluria gene panels.
- Diagnostic findings for Primary Hyperoxaluria panel were all due to homozygous or compound heterozygous pathogenic variants in the AGXT gene.
- Diagnostic findings for the nephrolithiasis panels:
 - The majority were identified with biallelic pathogenic variants in autosomal recessive genes: *GRHPR*, *ATP6VOA4*, *SLC7A9*, *CLDN16* and *ATP6V1B1*
 - One X-linked gene *CLCN5*
 - One autosomal dominant gene *CASR*
- All disease-causing variants detected have been reported in literature previously, supporting that genetic underpinnings play an important role in the etiology of kidney stones.
- Potential diagnostic findings:
 - An additional 6% cases (7 total) with variant(s) of uncertain significance that may potentially become diagnostic based on inheritance. Among these, half are autosomal recessive and half are autosomal dominant.
 - There are 2% cases (2 total) with pathogenic/likely pathogenic variant(s) in genes that have been reported as both autosomal dominant and autosomal recessive inheritance.
 - Familial segregation studies would be the next step to further investigate the potential pathogenicity of these variants and/or diagnosis of disease.

Methods

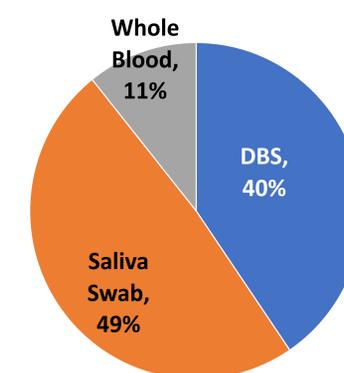
- In total 121 samples between June 2020 and February 2022 were tested at our laboratory.
- Sample types includes dried blood spot (DBS), saliva swab, and whole blood
- Samples came from three different types of clinics
 - Urology
 - Nephrology
 - Pediatric nephrology
- There are 55% males and 45% females.
- Next generation sequencing and deletion/duplication analysis were performed using a capture-based library on Illumina sequencer.

Client practice category



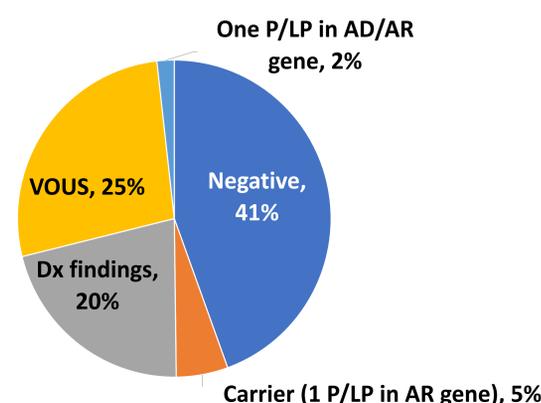
■ Urology ■ Pediatric Nephrology ■ Nephrology

Sample type distribution



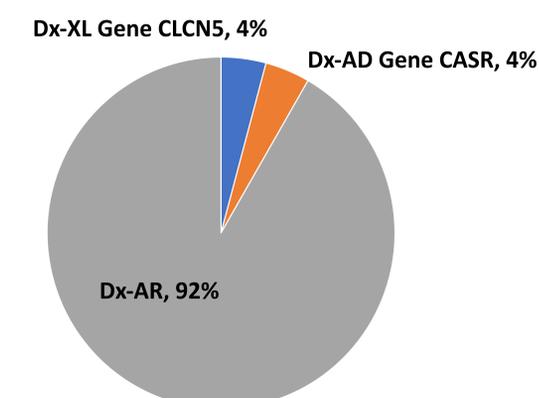
■ DBS ■ Saliva Swab ■ Whole Blood

Sequencing results/variant types



■ Negative ■ Carrier (1 P/LP in AR gene)
■ Dx findings ■ VOUS
■ One P/LP in AD/AR gene

Inheritance of diagnostic findings



■ Dx-XL Gene CLCN5 ■ Dx-AD Gene CASR ■ Dx-AR

CONCLUSION

- Our 35-gene Nephrolithiasis and 3-gene Primary Hyperoxaluria sequencing panels have a high diagnostic yield with significant potential for clinical use.
- Given the prevalence of genetic causes in these diseases, gene sequencing should be considered as first line testing and standard of care.
- Familial segregation study is the next step to reach possible diagnosis for those identified with variant(s) of uncertain significance or one pathogenic/likely pathogenic variant with uncertain inheritance.

Scan to Download

