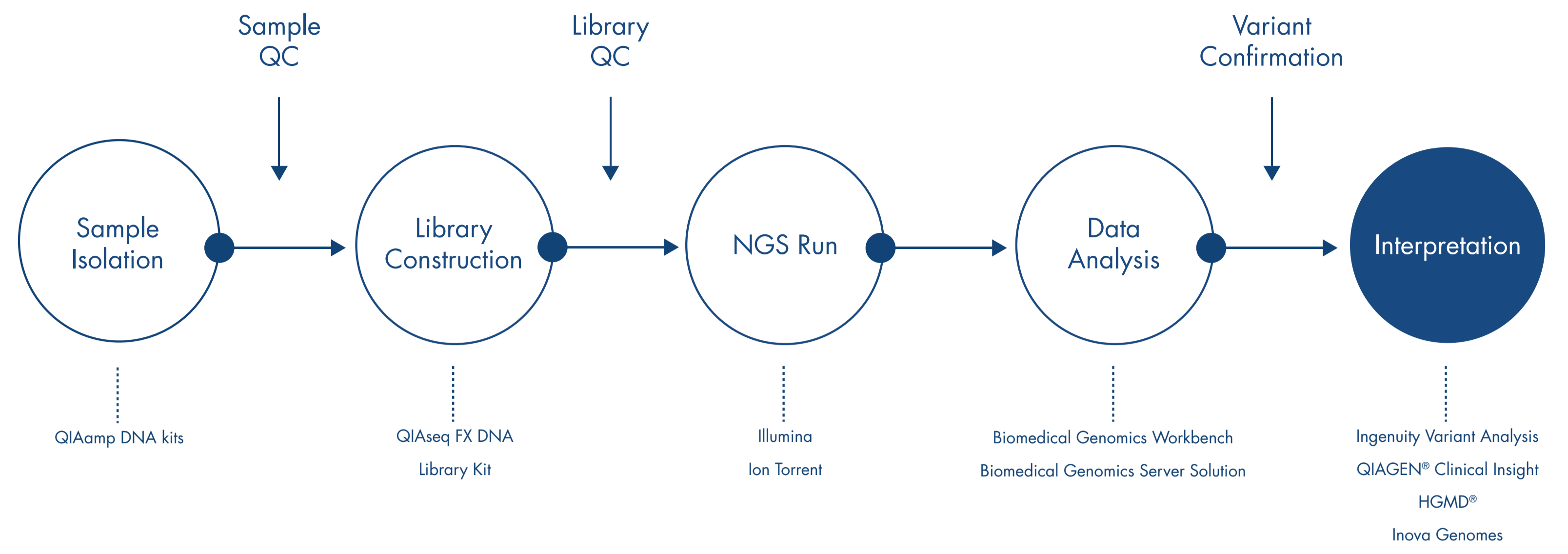
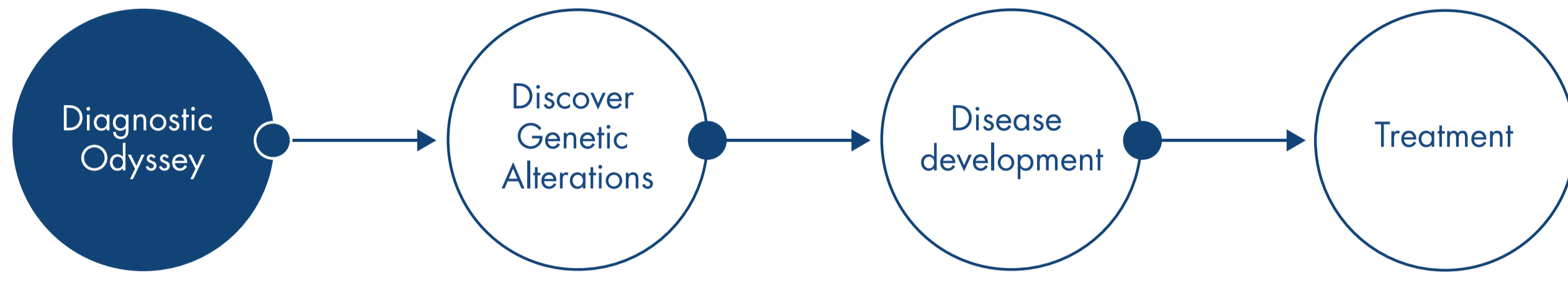
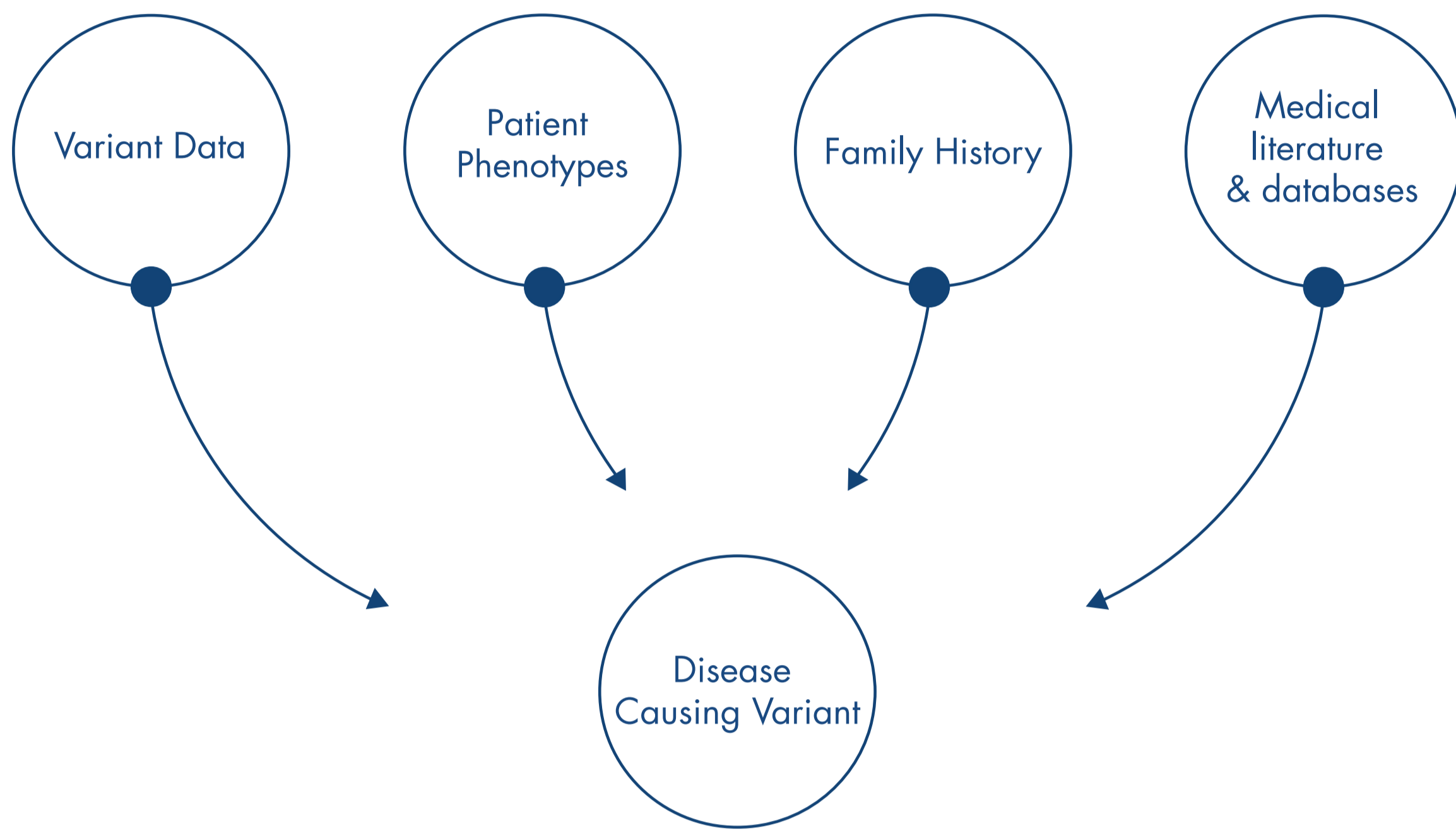


Clinical Genome and Exome Sequencing (CGES)



Identify Disease Causing Variants



Causal Variant Discovery & Interpretation

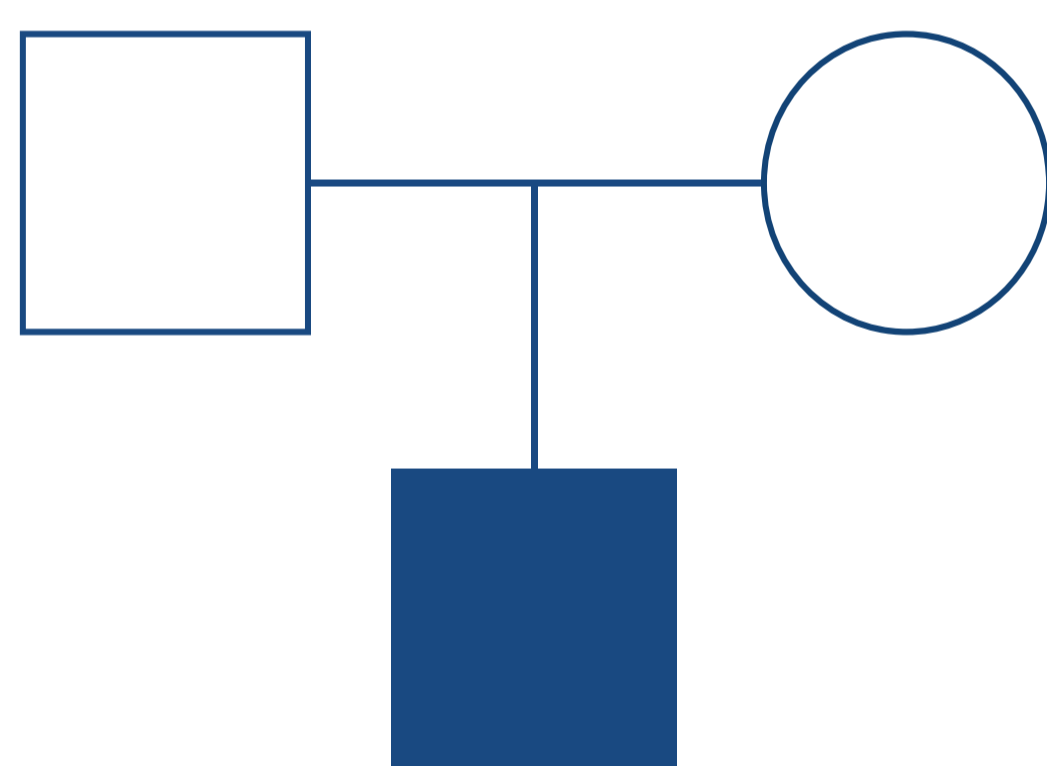
Confidence Call Quality > 20
 Variants not present in highly variable exonic regions

Common Variants MAF < 0.5
 1000G, NHLBI-EVS, ExAC, Allele Frequency Community

Predicted Deleterious Pathogenic or Likely Pathogenic (ACMG)
 HGMD
 LoF and Missense

Genetic Analysis Variant shared by proband and mother

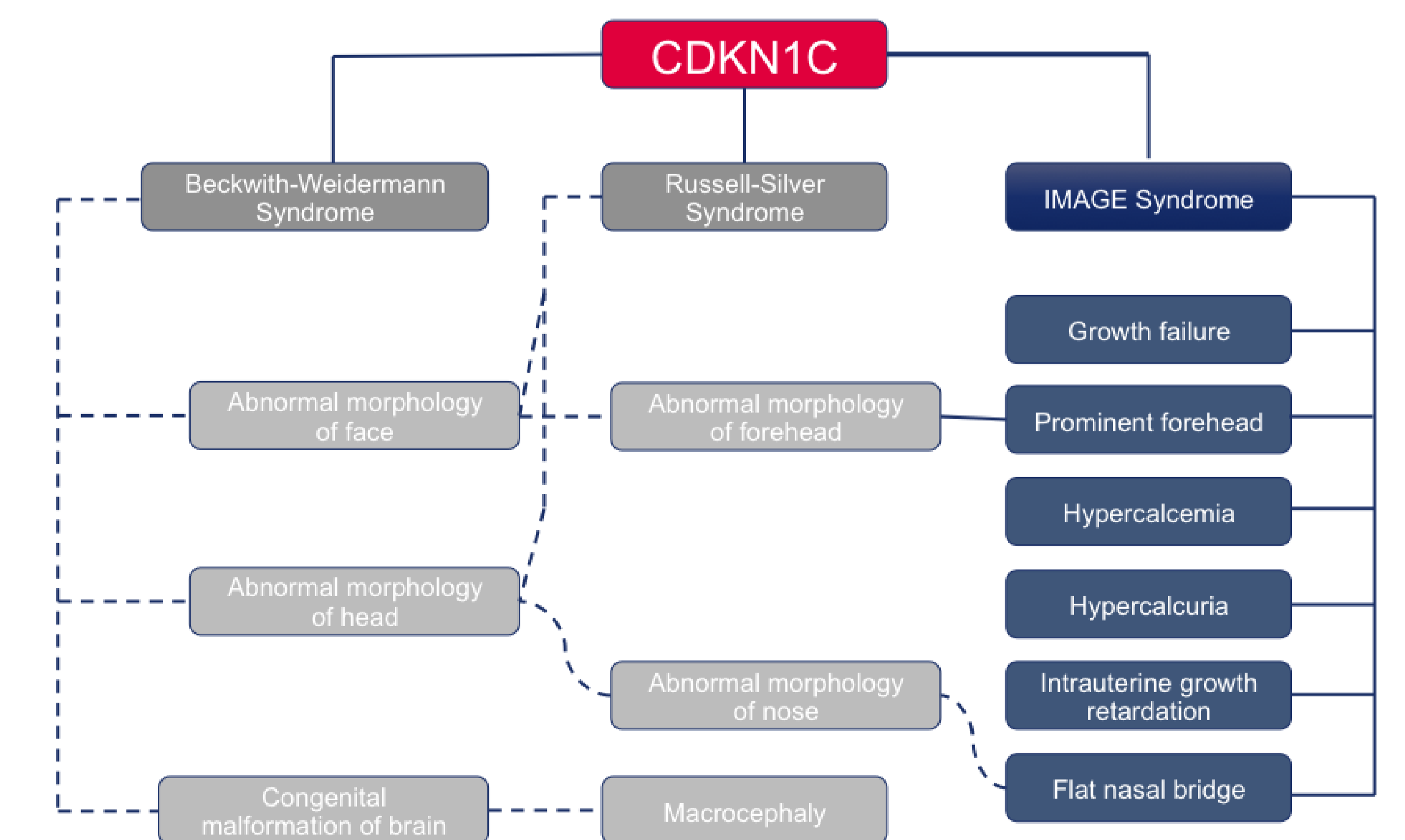
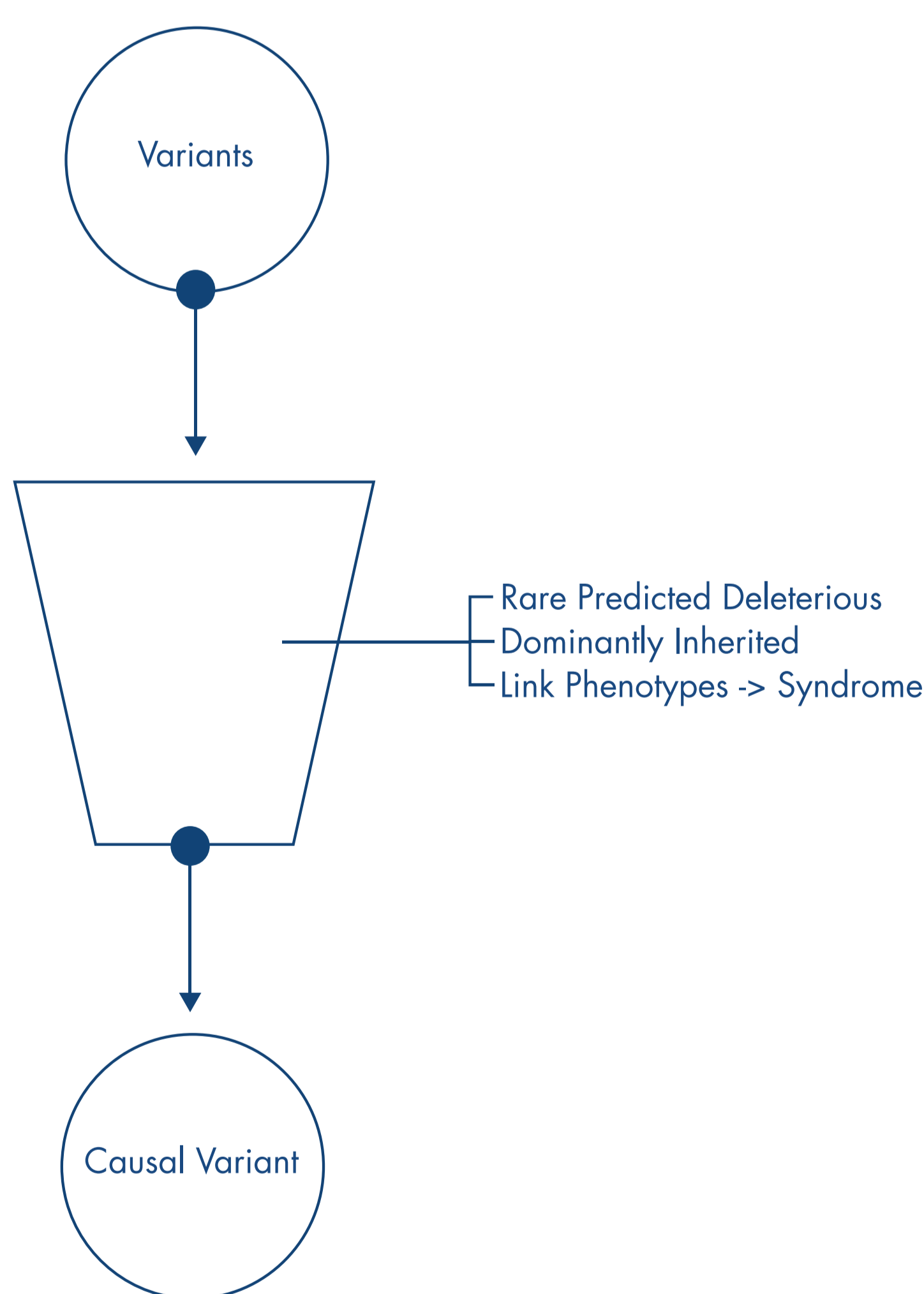
Biological Context Phenotypes-
 Prominent forehead; Macrocephaly; Hypospadias; Intrauterine growth retardation; Flat nasal bridge



- Phenotypes**
- Prominent forehead
 - Macrocephaly
 - Hypospadias
 - Intrauterine growth retardation
 - Epiphyseal dysplasia
 - Metaphyseal dysplasia
 - Flat nasal bridge

Syndrome Inference

Disease	Gene	Chr	Position	Causal	MOI	Genotype	Phenotype
IMAGE syndrome	CDKN1C	11	2905353	yes			
Donnai-Barrow syndrome	LRP2			yes	AR		
Smith-Lemli-Opitz syndrome	LDLR			no	AR		
Leprechaunism	INSR			yes	AR		
Seckel syndrome 1	CENPJ			yes	AR		
MELAS syndrome	MT-ND1			yes			
Lennox-Gastaut syndrome	MT-ND1			no			



Summary

- Fastest & most accurate application with built-in false positive check
- Leverage peer-reviewed literature content and network analysis to infer syndromes from patient phenotypes
- Discover known and novel causal variants