

Precision medicine, for sharper decisions.

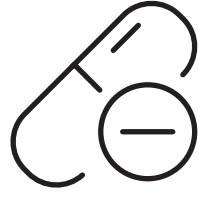
Why precision medicine is superior



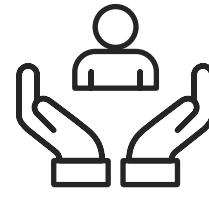
Better insights to disease drivers



Improved safety and efficacy profiles



Avoided costly drug attrition



Tailored therapies for the best outcomes

Traditional approaches



One size fits all



25% of cancer patients respond¹

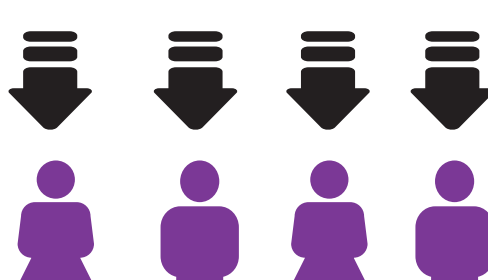
Precision medicine



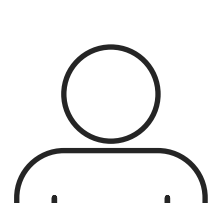
Stratification



Tailored therapy



100% patient response



The right individual



The right drug

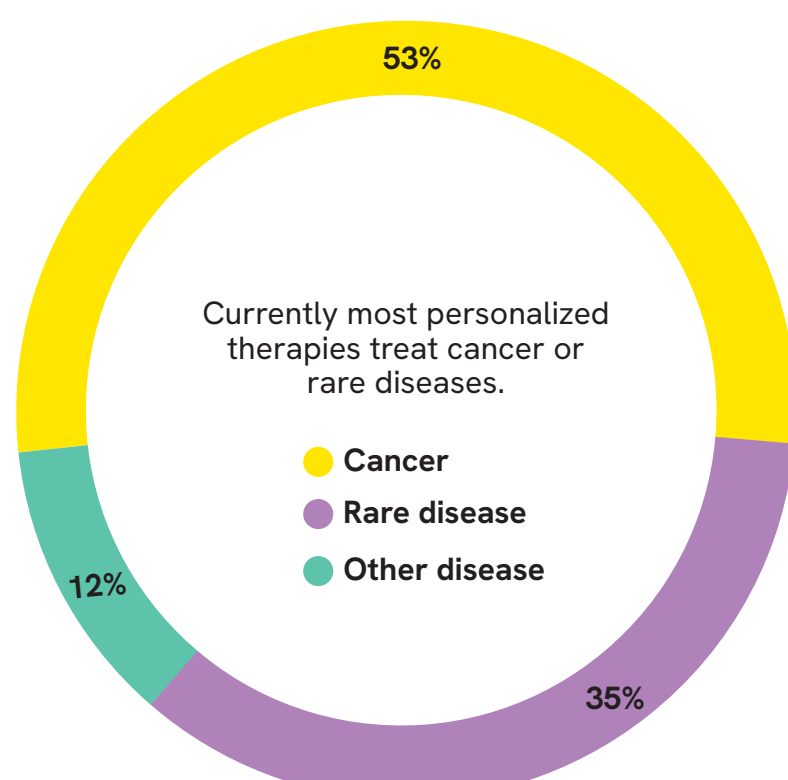
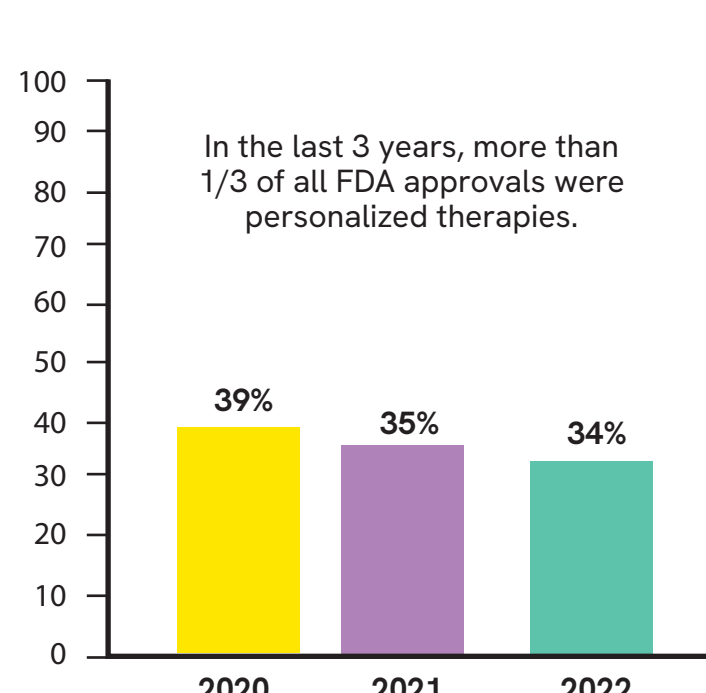


The right dose

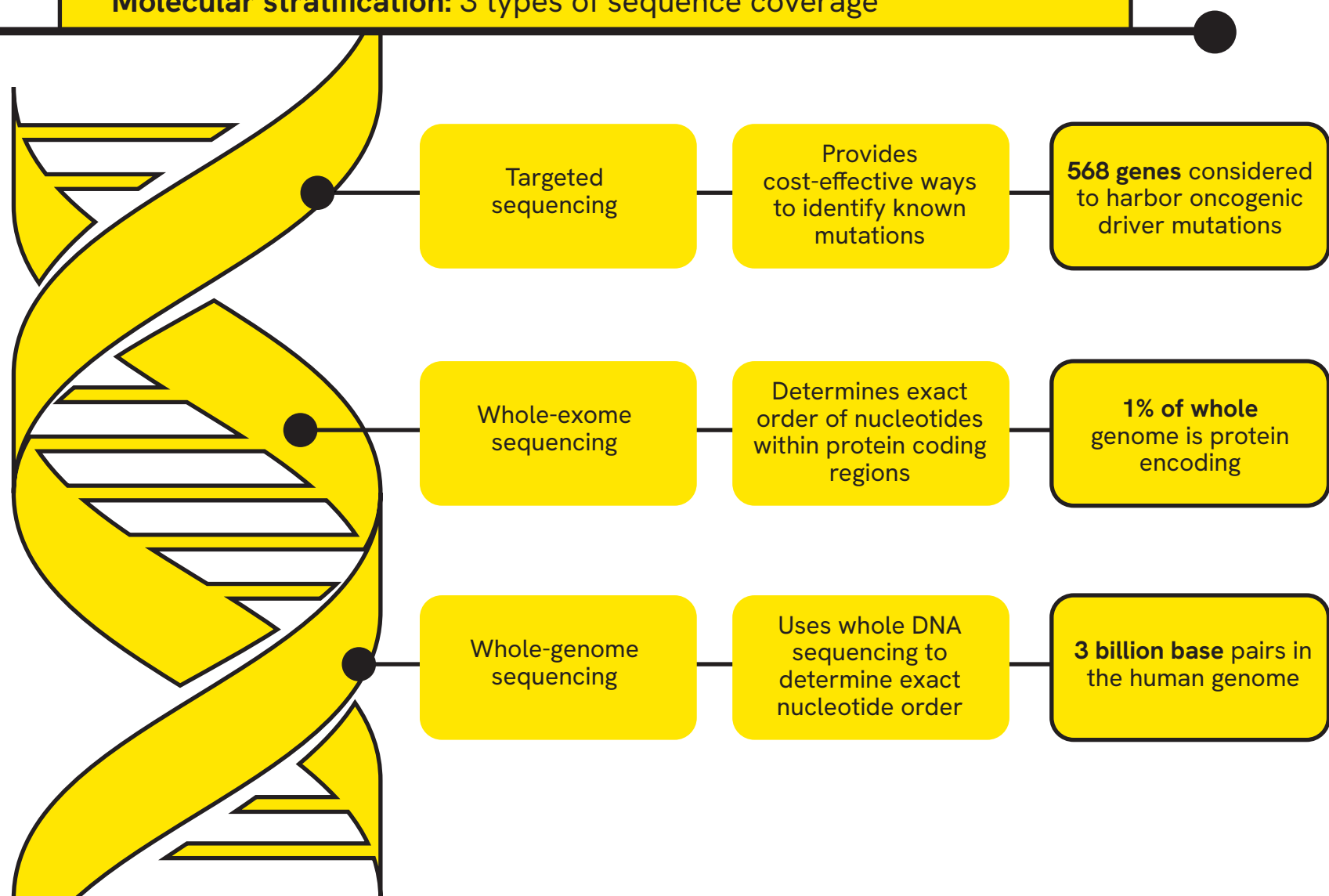


The right time

Clinical success with precision medicine²



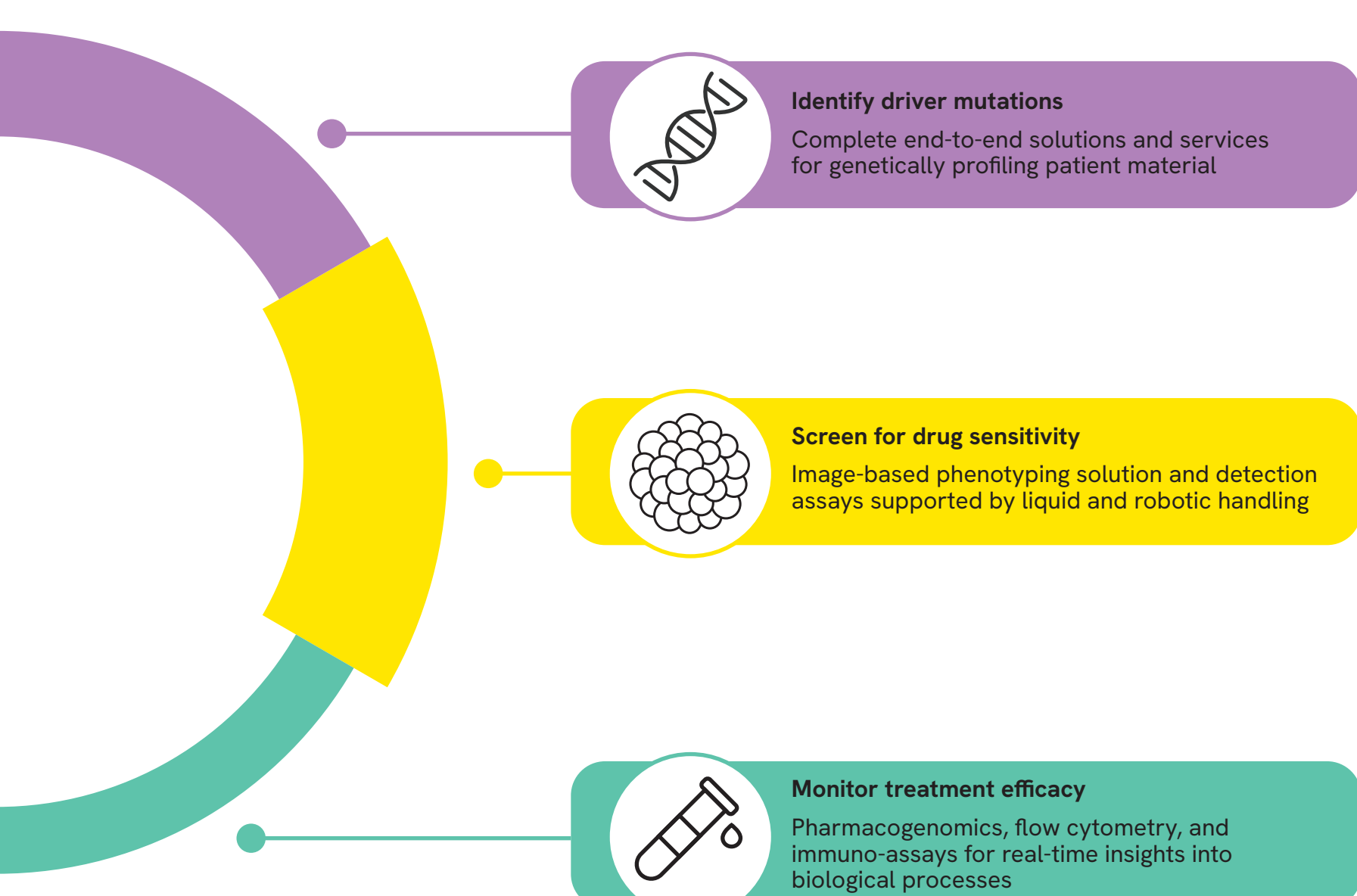
Molecular stratification: 3 types of sequence coverage



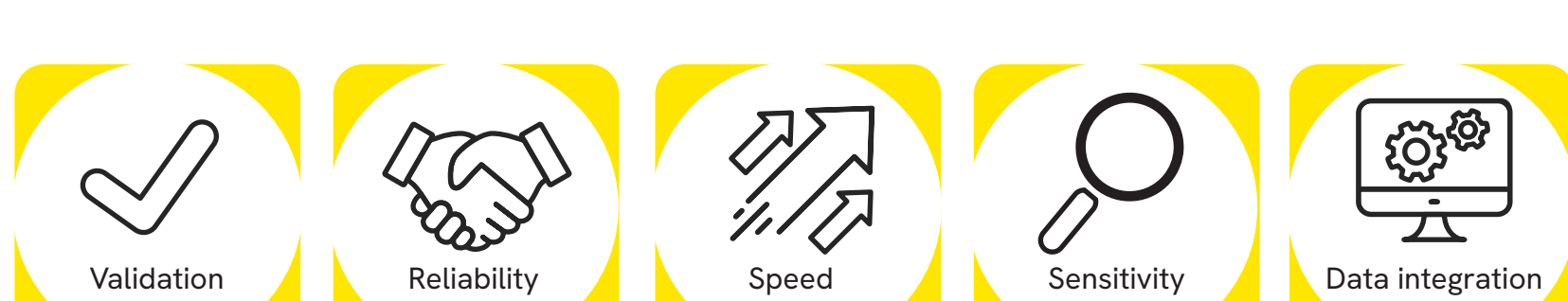
Functional stratification: which ex vivo model system is right for you?

| | Direct profiling | Cultured models | PDX models |
|-------------|--|---|--|
| Pros | Rapid assessment post excision retains tumor expression and immune populations | Expansion of primary tissue in physiologically relevant format enables great number of drug combinations to be tested | Replicates complex <i>in vivo</i> tumor stromal environments |
| Cons | Limited material means few drug combinations can be tested | Difficult to create and introduces risk of not faithfully recapitulating patient material | Results only available after months of culture; question over ethical responsibility |

Using knowledge to build precision



Achieving stratification clarity



Find the right tools for success at www.revvity.com/precisionmedicine

¹<https://hitconsultant.net/2014/04/03/infographic-the-rise-of-personalized-medicine/>

²The Scope & Significance of Progress in 2022 Personalized Medicine at the FDA