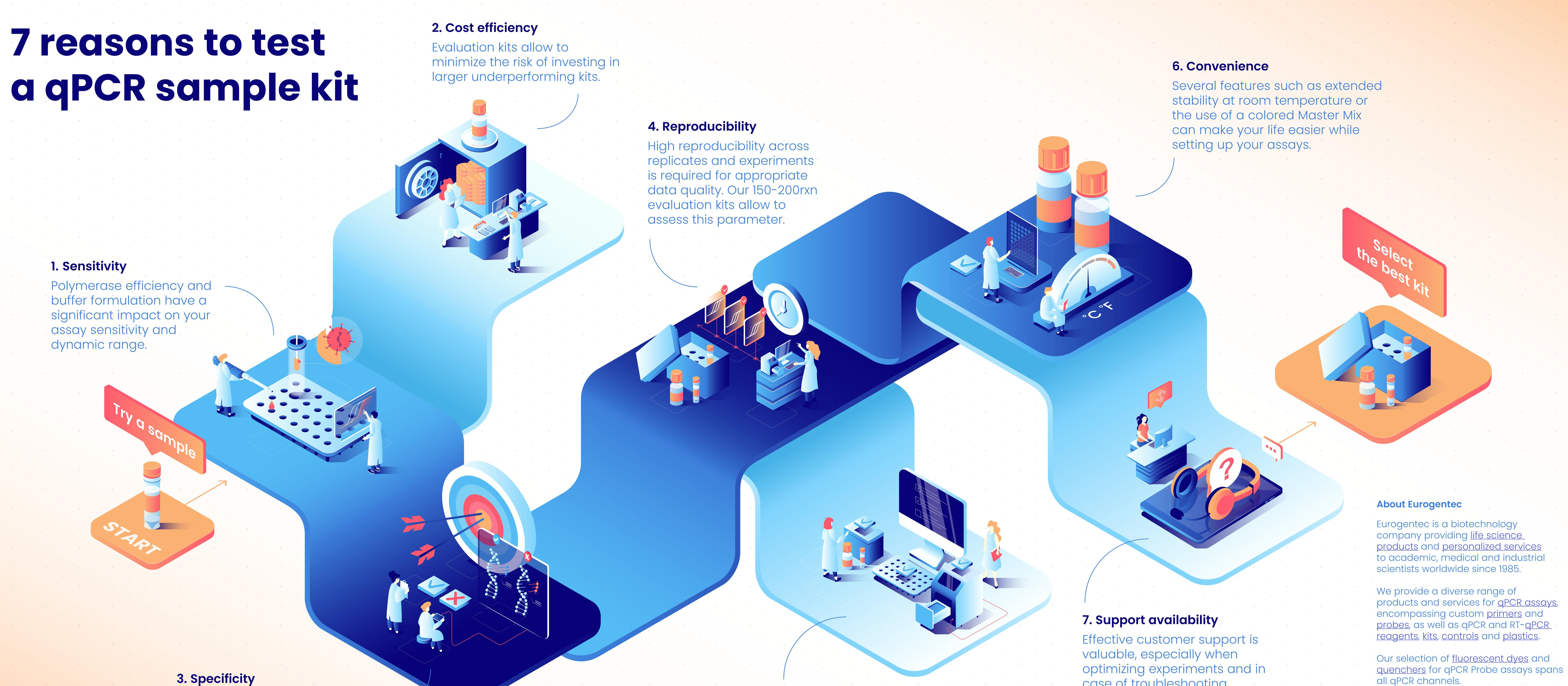


Some qPCR kits may require optimization for some targets or

applications while others don't.



5. Compatibility

Ensures that the kit is

samples, assays, and

compatible with your specific

instrument (e.g. ROX level).

In addition to catalog items, we produce <u>custom amplification reagents</u> and offer <u>custom kit assembly</u> services.

reproducible results.

When paired with our <u>Takyon® qPCR</u> or

RT-qPCR Master mixes, they ensure the

achievement of sensitive, accurate, and

case of troubleshooting.