



Identify targets others miss.

Pooled CRISPR screening solutions, from
basic research to lead optimization.

Leverage pooled screening to simultaneously target multiple genes within a cell population. Access solutions designed to help you reduce costs while maintaining the scalability and efficiency required for modern functional genomics research.

We're here to empower you with reliable reagents and robust platform solutions tailored for pooled screening workflows. Our end-to-end systems enable high-throughput research with exceptional precision, accelerating the identification and validation of novel drug targets in key areas such as oncology, neurodegeneration, and rare diseases.

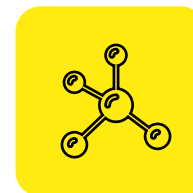
Whether you're identifying essential genes, uncovering gene interactions, or mapping complex biological pathways, our integrated solutions streamline your journey from target discovery to therapeutic development—bringing innovative treatments to patients faster.



Basic
research



Target
discovery



Target
validation



Lead
identification



Lead
optimization

