



# Your pathway toward CRISPR workflows.

Arrayed CRISPR screening,  
a better understanding starts here.

Gain a clear phenotype to genotype correlation with solutions to help perform CRISPR-based perturbation within separate wells of a microplate. Identify and validate hits and understand off-target effects of your drug candidates.

Providing solutions to automate high-throughput workflows, help maintain quality control, integrate large-scale genomic data for meaningful insights, and bring cost-effectiveness to your processes.

Improve your arrayed screening capabilities and gain the advantages of:

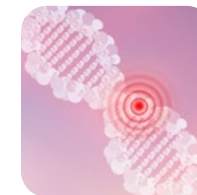
- Multi-parametric read-outs for cell viability, morphology, differentiation, and proteomic profiling
- Ability to interrogate more complex cell models
- Greater choices in reagent formats, including non-viral transfection and cherry-picking options
- Clear phenotype to genotype correlation
- Improved editing efficiency by targeting the same gene with several gRNAs in one well



Disease  
models



Pathway  
analysis



Target  
identification

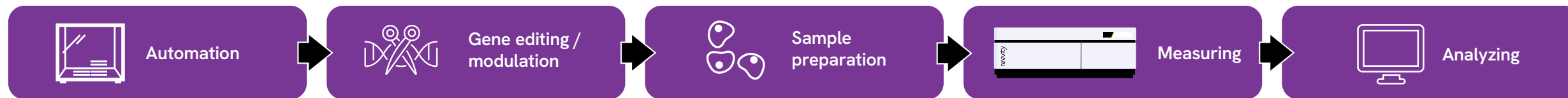


Target  
validation



Cell and gene  
therapy

## Workflow



## Revvity solutions

### Automated liquid handling workstations

Minimize errors, reduce hands-on time, and increase throughput and reproducibility.

### Dharmacon™ reagents

Gene editing reagents that offer flexibility to achieve any edit, in any cell type, at any scale.

### PhenoVue™ cellular imaging reagents

Optimized & verified reagents including cell painting kits, organelle and cell compartment stains, and cell function reagents.

### Multimode microplate readers

Plate readers equipped with the latest detection technologies.

### Cellular imaging software

Acquire, visualize, analyze, and discover with the Harmony™ imaging and analysis software and the Signals Image Artist™.

### BioLegend Zombie fixable viability kit

Stain with non-permeant dye for live cells but permeant to cells with compromised membranes.

### Engineered cell lines

Ready-to-go validated knockout models, point mutations in cancer-relevant cell lines, and Cas9 and dCas9-VPR cell lines.

### Immunoassays

Offering rapid, homogeneous, and easy to use AlphaLISA™ and HTRF™, no wash assays.

### PhenoPlate microplates

Optimized for high-content imaging and engineered to deliver superior images and high-quality data.

### Revvity Signals

First unified, cloud-native SaaS R&D solution that streamlines scientific complexity and promotes collaboration.

### Cell counters

High-throughput cell counting that uses a multi-well microplate-based format.

### BioLegend LEGENDplex™

Quantitate up to 14 analytes with flow cytometry-based multiplex immunoassays.

### High-content imagers

Opera Phenix™ Plus and Operetta™ CLS high-content imaging systems feature confocal spinning disk technology for exceptional image quality.

### Functional genomics screening services

Arrayed screens using CRISPR, RNAi, and Pin-point™ base editing technologies.

### MojoSort™ magnetic cell sorting

Isolate and purify cells from heterogeneous populations.

revvity