

# Complement your screening with scalable transcriptomics.

## The challenge:

Phenotypic screens show what happened, but not always why.

### MoA blind spot

Phenotypic readouts rank hits by severity but often provide limited insight into mechanism

### Hit prioritization cap

Without transcriptional context, morphological hits advance with similar confidence, increasing risk of pursuing off-target or low-value compounds.

### Off-target risk undetected

Imaging and viability may not capture pathway-level off-target activity. Gene-expression signatures reveal unintended biology before *in vivo* studies.

## The solution:

Alithea's MERCURIUS™ DRUG-seq & Total DRUG-seq  
Built for scale.

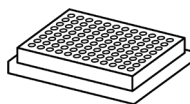
Revvity is an authorized distributor of:

ALITHEA  
GENOMICS



### Extraction-free

Direct from cells.  
Eliminates the need for RNA isolation.



### Early multiplexing

Sample barcoding enables pooling of up to 1,000+ conditions in a single tube.



### Genome-wide coverage

Gene-expression signatures across the genome.



### Ultra low-cost

Significantly cheaper than conventional alternatives.

## Two use cases in screening workflows

### Compound screening

- **MoA assignment:** cluster compounds by transcriptional signature to support MoA inference
- **Hit prioritization:** rank hits by pathway activity alongside phenotypic severity
- **Off-target insight:** flag potential unintended pathway activity before *in vivo* studies

**Morphology + transcriptomics =  
higher-confidence MoA, smarter hit triage.**

### CRISPR & siRNA screens

- 384-well arrayed CRISPR/siRNA screens with transcriptional readout.
- On-target confirmation: assess expected pathway-level transcriptional responses
- Off-target identification: detect potential unintended pathway activation or silencing
- Pair with imaging for dual phenotypic plus transcriptional annotation of every perturbation

**From genetic perturbation to pathway hypothesis  
in a single workflow**

## Selection guide

Available in 96-, 384-, and 1,536-well formats to match screening workflows.

	DRUG-seq	Total DRUG-seq
Data output	Gene-level expression signatures	Full transcriptome resolution
RNA captured	Poly(A) RNA (mRNA only)	Total RNA (mRNA + non-coding)
Transcript coverage	3' gene-level counting	Full-length transcript coverage
Reads/sample	~1-2M	~2-5M

- **DRUG-seq** enables high-throughput profiling with efficient gene-expression signatures for large-scale screens.
- **Total DRUG-seq** extends this approach to full transcriptome resolution, capturing isoforms and non-coding RNA for deeper biological insight.

## Did you know?

Revvity provides an end-to-end screening workflow of instruments and reagents to support MoA analysis, hit prioritization, and off-target assessment.

### See beyond the phenotype

Contact your Revvity representative to integrate DRUG-seq into your screening workflow.



revvity