

HIGH-THROUGHPUT SCREENING & LEAD OPTIMIZATION PRODUCTS

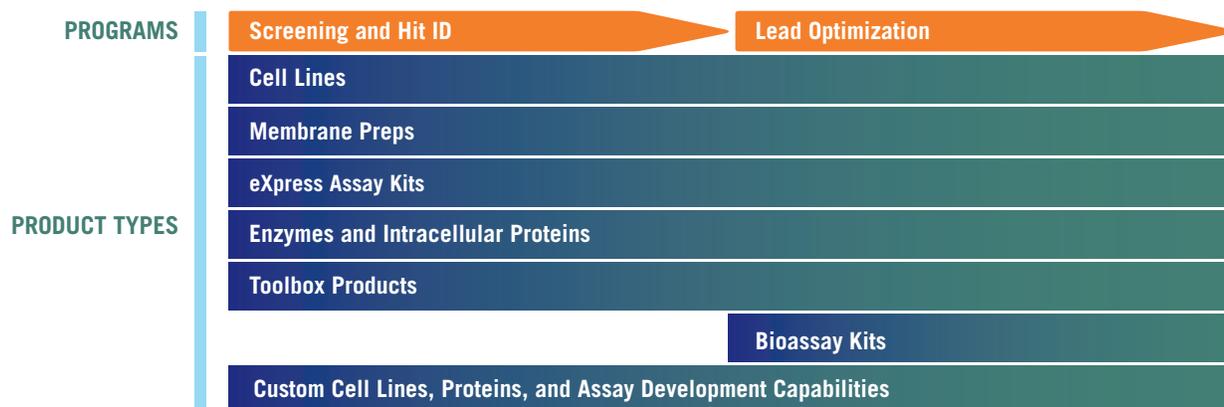
Robust Cell-Based Assays to Identify Therapeutic Hits Rapidly

SCALABLE, SIMPLE CELL-BASED ASSAYS FOR SCREENING HITS TO LEAD OPTIMIZATION

Developing therapeutics, small molecule or biologics, presents unique challenges at each development phase. Utilize Eurofins DiscoverX high-quality cell-based assays to support drug discovery and development programs from high-throughput screening through lead optimization

and beyond. Access the largest available portfolio of products including cell-based functional and binding assays, cell lines, and enzymes covering the top therapeutic target classes including GPCRs, Kinases, NHRs, Checkpoints, Ion Channels, and Cytokine Receptors.

Expedite your candidate through the drug discovery pipeline with Eurofins DiscoverX assays



SUPPORTING DRUG DISCOVERY & DEVELOPMENT PROGRAMS

- **Enabling Technologies** – Improve productivity and effectiveness of your screening, lead optimization, and SAR campaigns to accelerate the advancement of new therapeutics
- **Simple, Fast, and Scalable** – Rapid, homogeneous, and simple mix-and-read assay format that is scalable from 96- to 3456-well and does not require washing, centrifugation, or filtration
- **Robust Assays** – Cell-based assays with large signal-to-background ratio, high precision, and lot-to-lot reproducibility
- **Qualified and Validated** – Extensively optimized for hundreds of targets used for screening in billions of data points and cited in thousands of peer-reviewed publications



Visit discoverx.com to start successfully screening today!