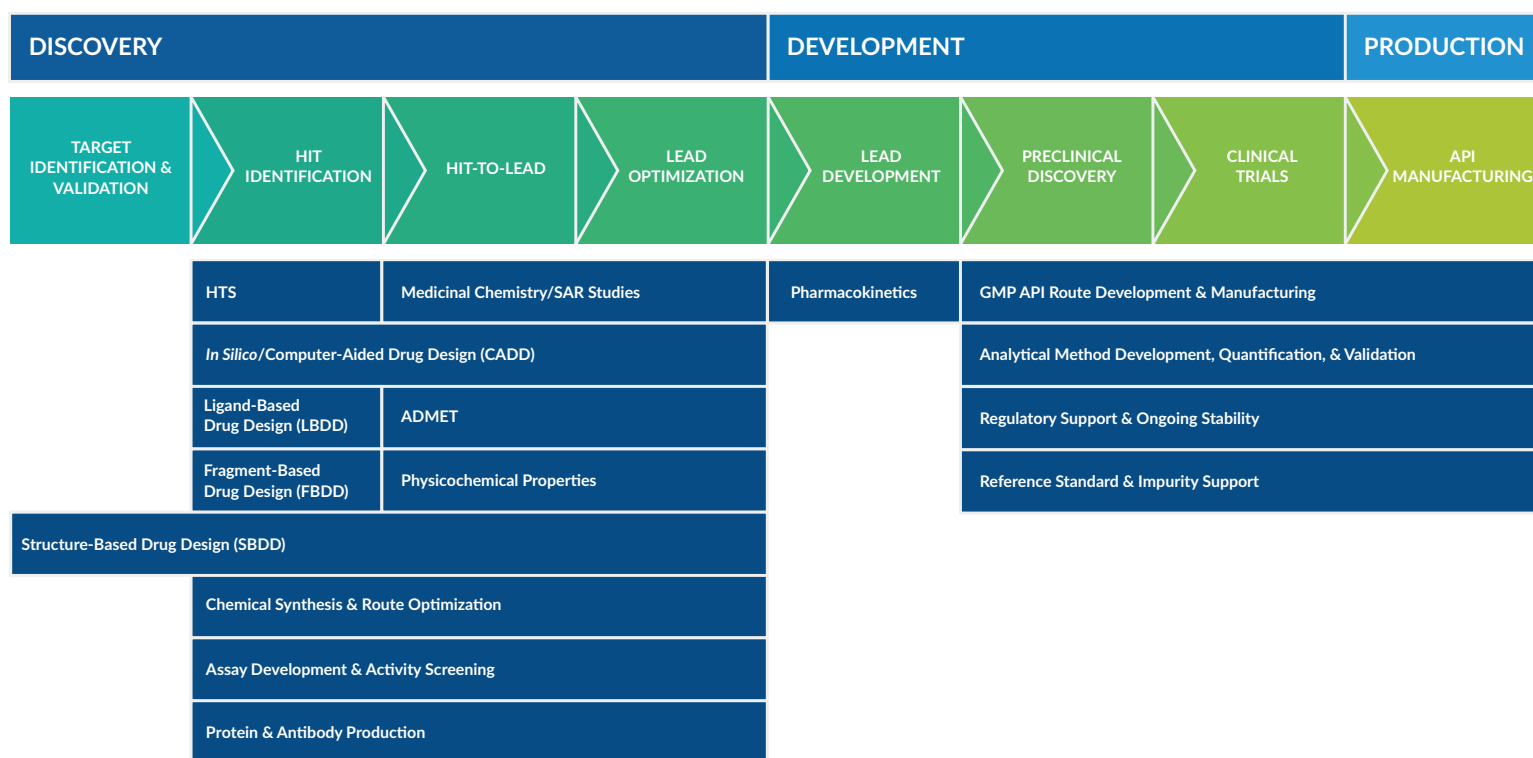


Integrated Drug Discovery

With our wide range of chemistry and biology services, Cayman is perfectly positioned to provide integrated, multidisciplinary development and preclinical drug discovery services to accelerate your journey from target to clinic. Our drug discovery services are powered by expertise in computational chemistry, compound screening, medicinal chemistry, chemical synthesis, structural biology, and GMP API manufacturing. Additional experts in assay development, cellular metabolism, immunology, and cell and molecular biology can develop custom assays and model systems to test your compounds in the most relevant contexts.

Drug Discovery & GMP API Manufacturing Services



Why Partner with Cayman?



Client-centric Approach

- Your needs are our top priority



Scientific Expertise

- Skilled across drug & target classes



State-of-the-art Facilities

- In-house platforms & trusted partners



Comprehensive

- Broad portfolio of products & services



Flexible

- Single-point or end-to-end support



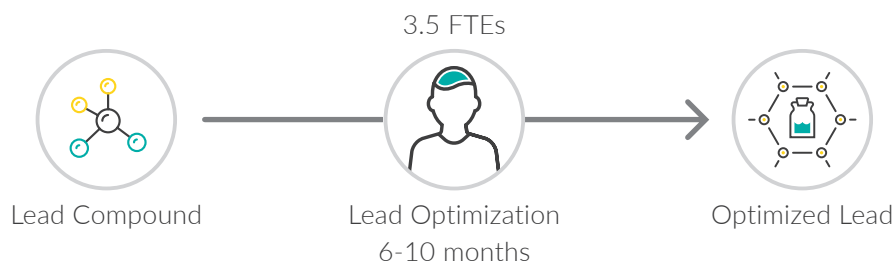
Quality

- Validation & controls at every step

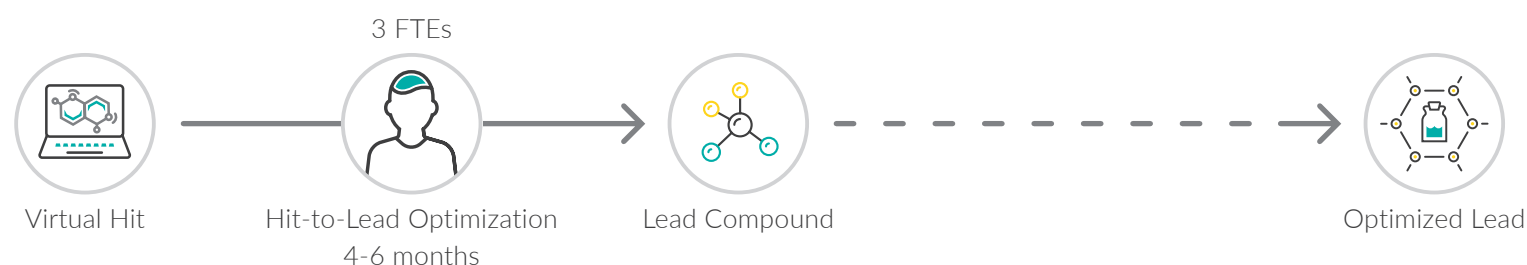
Our Approach

We challenge the lengthy, costly, and laborious process of building and testing thousands of compounds by adopting an intelligent compound design and selection approach guided by the expertise of our team. We emphasize quality over quantity. Each compound is designed to answer a structure-activity relationship (SAR) question to minimize the synthesis of fruitless compounds. We utilize a variety of computational, structural, and biophysical methods in our state-of-the-art facilities to accelerate and streamline drug design and discovery.

Lead Series



Backup Lead Series



Learn how we accelerated the discovery and design of novel heterocyclic inhibitors from hit identification through lead optimization in less than 10 months.

Download the poster at www.caymanchem.com/heterocycleinhibitors

Medicinal & Computational Chemistry

- Hit identification with virtual high-throughput screening (vHTS)
- Hit-to-lead optimization with iterative SAR studies
- Activity screening, selectivity, & potency

Structure-based Drug Design

- *In silico* drug design/computer-aided drug design (CADD)
- Structure determination by X-ray crystallography and cryo-EM
- Surface plasmon resonance (SPR) with Biacore™ 8K, thermal shift assay (TSA), and static or dynamic light scattering (SLS/DLS)

Chemical Synthesis

- Small molecules, drug-like heterocyclic compounds
- Physicochemical property testing
- Lipid nanoparticle development

GMP API Manufacturing

- Process scaling, route optimization, and pre-GMP process development
- Preclinical/clinical API synthesis and method validation
- US-based CGMP API manufacturing

Let's discuss your project needs at www.caymanchem.com/services