

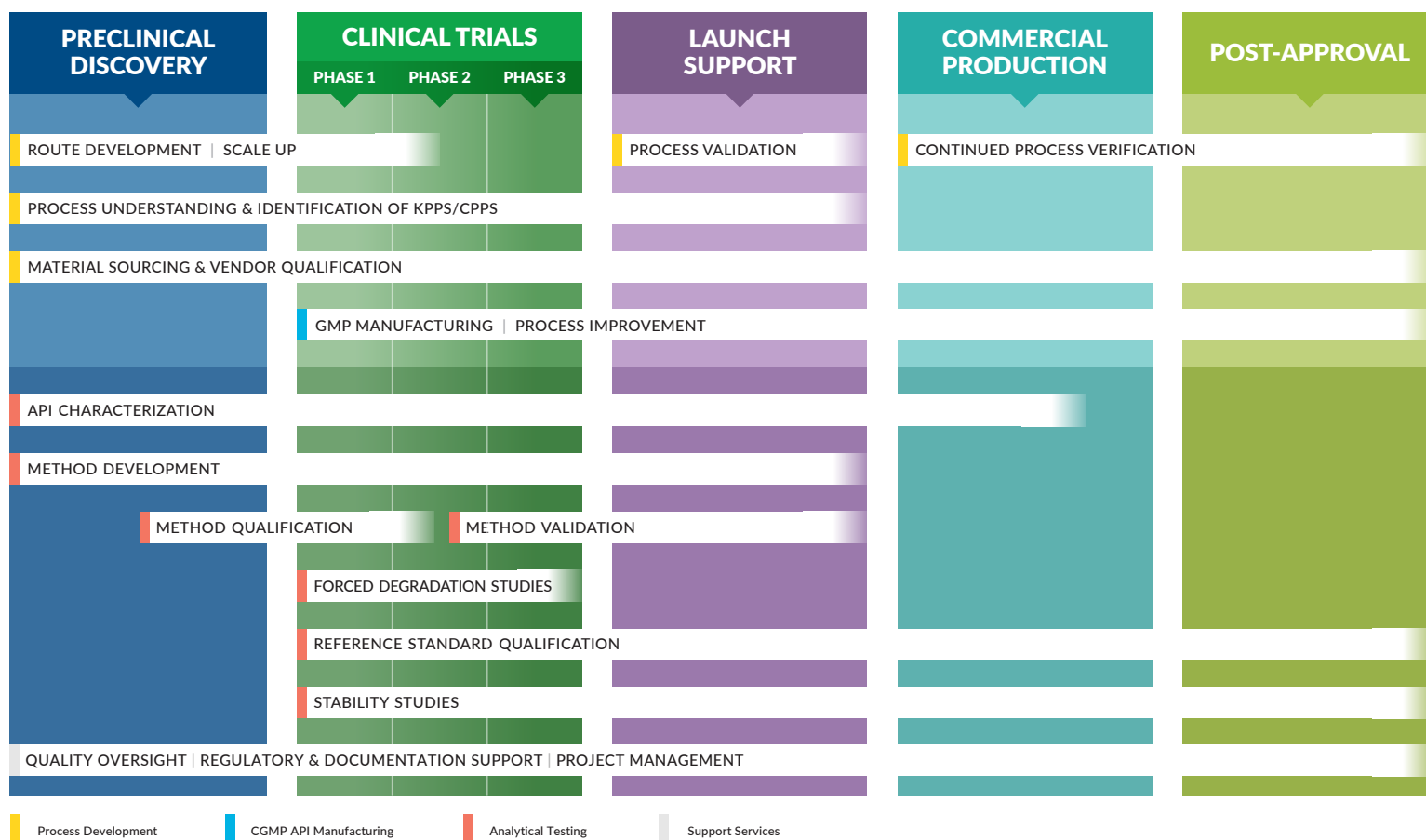
CGMP Services

A Trusted Partner from Molecule to Market

Cayman is a global contract development and manufacturing organization (CDMO) that supports academic, pharma, and biotech clients across the entire drug development lifecycle, from drug discovery through clinical trials to commercialization. With over 40 years of experience in API manufacturing, we have proven expertise in API development and manufacturing in full compliance with CGMP standards. We offer a collaborative, flexible approach, allowing our clients to create a customized service plan that meets their needs.



Our Services Include:



Chemistry Capabilities

- Prostaglandins & Prostanoids
- Nucleoside Analogs
- Aminoisoquinoline & Aminoisoquinolone Analogs
- Tryptamines & Other Psychedelics
- Fixed & Ionizable Cationic Lipids
- Heterocyclic Compounds
- Custom Molecules Upon Request

Cayman is proud to be one of the first CDMOs offering psychedelics for use in human clinical trials

20 years of experience in navigating the regulatory landscape of manufacturing, importing, and exporting APIs and DEA Schedule I-V controlled substances.

Ability to work with formulation CDMOs and the DEA to secure quota and deliver psilocybin, psilocin, and other psychedelic APIs.

Our Excellence

Process Development

Our highly skilled chemists specialize in fit-for-purpose route design of complex, multi-step organic syntheses. They leverage the extensive R&D capabilities and diverse chemistry experience of Cayman Chemical to improve process development and accelerate innovation.

CGMP API Manufacturing

Our expertise ensures a smooth transition during production scale-up that maintains process efficiency and consistent product purity. Our dedicated CGMP production space supports API production from multi-gram to multi-kilogram quantities, and we have a proven history of CGMP compliance during all phases of API manufacturing.

Analytical Testing

Cayman provides comprehensive analytical method development, qualification, and validation services to ensure product quality, consistency, and safety during all phases of API manufacturing. Our analytical team customizes phase-appropriate analytical testing programs to tailor to the individual requirements of each API while conforming to stringent regulatory requirements.

Support Services

We provide end-to-end support services at every stage to help advance your drug development program, including quality oversight, regulatory and documentation support, and a PMP-certified project manager who is your single point of contact for information and communication through the drug development cycle.

What Clients Are Saying

Cayman Chemical places paramount emphasis on meeting customer needs and fulfilling their specific requirements. The manufacturing site stand out for its impeccable cleanliness, efficient organization, and appropriate equipment setup, particularly tailored for the production of APIs.

Contract Auditor

Rephine