



Automating the Future of Synthetic Biology

Telesis Bio + Beckman Coulter Life Sciences

A Breakthrough Collaboration in DNA & mRNA Synthesis

Telesis Bio and Beckman Coulter Life Sciences have joined forces to revolutionize synthetic biology. By integrating the Gibson SOLA® Enzymatic Synthesis Platform with the fully automated Biomek Echo One Gene Synthesis Bundled Solution, powered by the Biomek i7 Hybrid Workstation and Echo 525 Liquid Handler, this collaboration delivers high-fidelity, scalable DNA and mRNA synthesis directly into your lab. The result is a streamlined, end-to-end workflow that accelerates innovation from design to synthesis with unmatched accuracy, reproducibility, and sustainability. Embrace non-toxic, aqueous chemistry and seamlessly incorporate third-party or custom constructs, all while gaining greater control and speed in your research without compromise.



Digital Sequence to Biological Endpoint in 1 Day. The Gibson SOLA® Platform on the Biomek Echo One System is a "digital-to-biological converter" that generates DNA or mRNA in less than one day.

Automated Workflow at a Glance

- Fully automated and requires only one user interaction for deck setup, offering complete walk-away capability.
- Divided into five modular method sections and includes safe stop points for flexibility.
- Workflow can be run as a full process or in discrete steps, depending on user needs.
- Entire synthesis process completes in under 24 hours, significantly faster than traditional service providers.
- SAMI Software supports scaling across up to four synthesis families with advanced scheduling and orchestration for high-throughput workflows.

The Gibson SOLA® Platform: A Digital-to-Biological Converter

This platform uses a hybridization-based, block assembly method that sets it apart from traditional phosphoramidite or TDT synthesis. It enables:

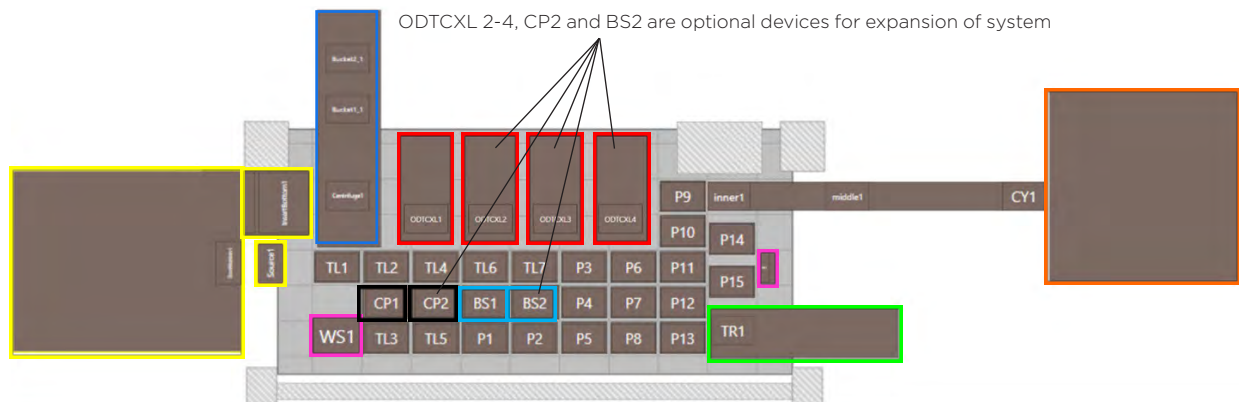
- Production of 10s to 100s of kilobases of high-quality DNA
- Over 80% full-length sequence yield
- Synthesis time reduced from weeks to under 24 hours

- Elimination of the need for dedicated HPLC columns
- On-demand, in-house synthesis of high-fidelity DNA with error rates <1:70,000 and no need for custom reagents

Biomek Echo One Gene Synthesis System: Automation That Delivers

This system integrates:

- Echo 525 Acoustic Liquid Handler, which enables non-contact, nanoliter-scale dispensing for rapid pooling and normalization
- Biomek i7 Hybrid Workstation, offering full walk-away capability with advanced pipetting, thermal cycling, and labware handling
- Biomek Software, providing intuitive setup, real-time monitoring, and error handling
- SAMI Software, supporting workflow scalability with advanced scheduling and orchestration
- Smart features such as Guided Labware Setup (GLS), DeckOptix Final Check (DFC), and Method Options Selector (MOS) for customizable workflows



The Biomek Echo One System deck for the Gibson SOLA® Platform that includes a total of four ODTC Thermocyclers.

Biomek i7 hybrid workstation with integrated Echo 525 acoustic liquid handler (yellow), ColdPlate (QInstruments, CP1 & CP2, black), BioShake D3000-elm (QInstruments, BS1 & BS2, blue), Automated Microplate centrifuge (Agilent, blue), on-deck thermocycler (Inheco, ODTCX1-4, red), Cytomat 2C (Thermo Fisher, orange), Biomek Wash Stations (pink), and High Capacity Trash (green).

About the Innovators

Telesis Bio

Pioneering synthetic biology with the Gibson SOLA® platform and BioXp® systems.

www.telesisbio.com

Beckman Coulter Life Sciences

A global leader in lab automation, empowering discovery since 1935.

www.beckman.com



© 2025 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries. ECHO is a trademark or registered trademark of Labcyte Inc. in the United States and other countries. Labcyte is a Beckman Coulter company. All other trademarks are the property of their respective owners.

For Beckman Coulter's worldwide office locations and phone numbers, please visit Contact Us at beckman.com

2025-GBL-EN-108620-v1

