

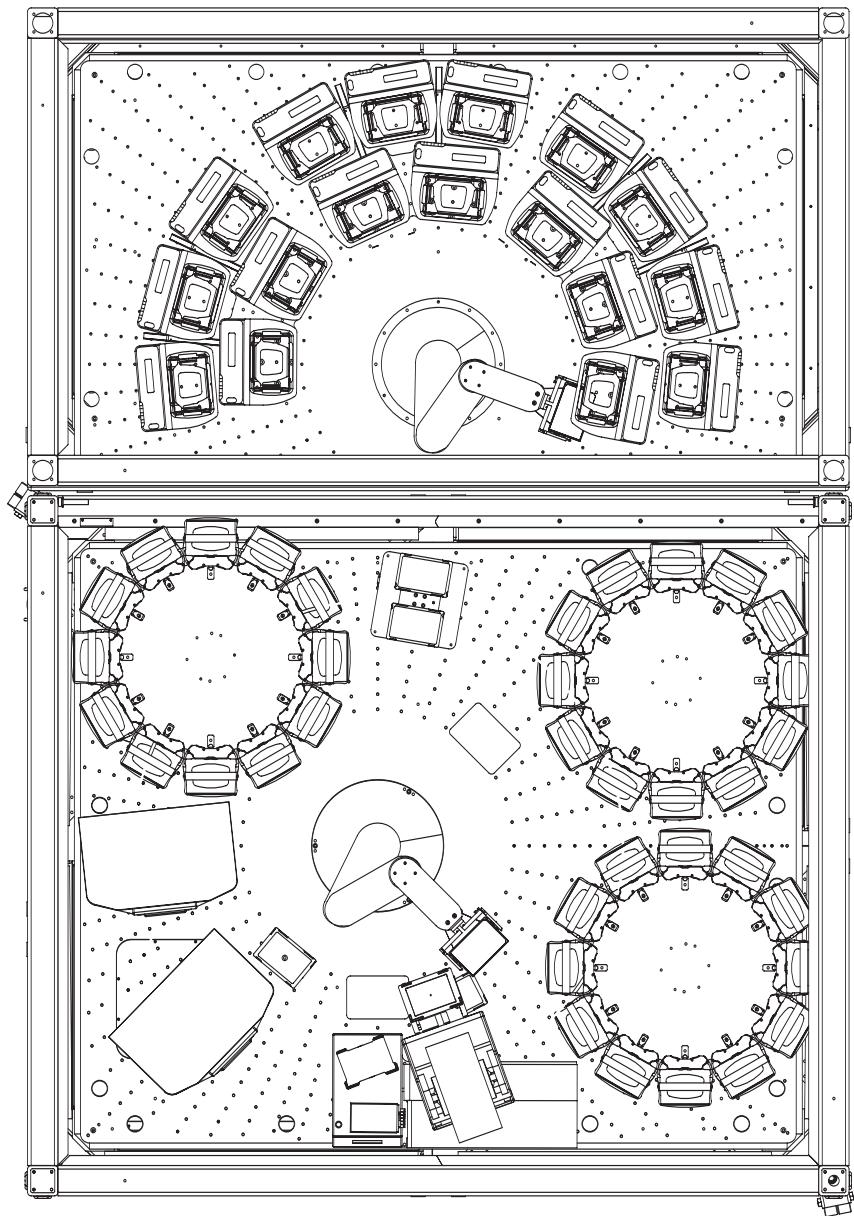
Agilent BioCel System Configuration

Compound Tube and Plate Replication

Application Bulletin

In compound management, the storage of compounds in individual tubes rather than in plates has great advantages — it allows individual access to compounds, easy reconfiguration of compound subsets and avoids capacity-consuming empty wells. This Agilent BioCel was designed for compound management to handle compounds stored in a variety of formats—Matrix 96 tube racks and Remp 384 microtube racks as well as standard microplates.

With this BioCel, the world of tubes (compounds) merges smoothly with the world of plates (assays). A high storage capacity for plates (17 Agilent Labware Stackers, 900 plates) and tube racks (3 random access carousels, 288 racks) allow the production of several sets of assay plates from compound tubes. The speed of the BioCel robotic arm allows the system to process its full plate capacity in less than 8 hours. The low volume, high precision pipettor Agilent Vertical Pipetting Station transfers volumes as low as 200 nL, aspirating and dispensing out of and into plates and tubes. Agilent VWorks scheduling software uses the Agilent Microplate Barcode Labeler to record and label barcodes on plates and tubes to ensure sample integrity. The BioCel can accommodate 1:4 and 4:1 tube/plate splits when moving between 96 and 384 (or 1536) formats. The Agilent PlateLoc Thermal Microplate Sealer seals the plates before they are stored in the output stackers. Tube sealing and/or environmental controlled storage for both tubes and plates could easily be accommodated into the system, making the system even more versatile.



Agilent Technologies

Module List

Component	Quantity	Function
Agilent robot	2	360° high speed robotic plate handler
Transfer station	2	Pass plates between robots
Agilent Vertical Pipetting Station	2	Low volume compound transfer
Agilent Microplate Barcode Labeler	1	Read and apply bar codes
Agilent PlateLoc	1	Thermal plate sealer
Carousel	3	Matrix tube rack access and storage
Agilent Labware Stacker	17	50-plate room temperature storage

www.agilent.com/lifesciences/automation

This item is intended for Research Use Only. Not for use in diagnostic procedures. Information, descriptions, and specifications in this publication are subject to change without notice.

Agilent Technologies shall not be liable for errors contained herein or for incidental or consequential damages in connection with the furnishing, performance, or use of this material.

© Agilent Technologies, Inc. 2009
Published in the U.S.A. February 25, 2009
5990-3622EN



Agilent Technologies