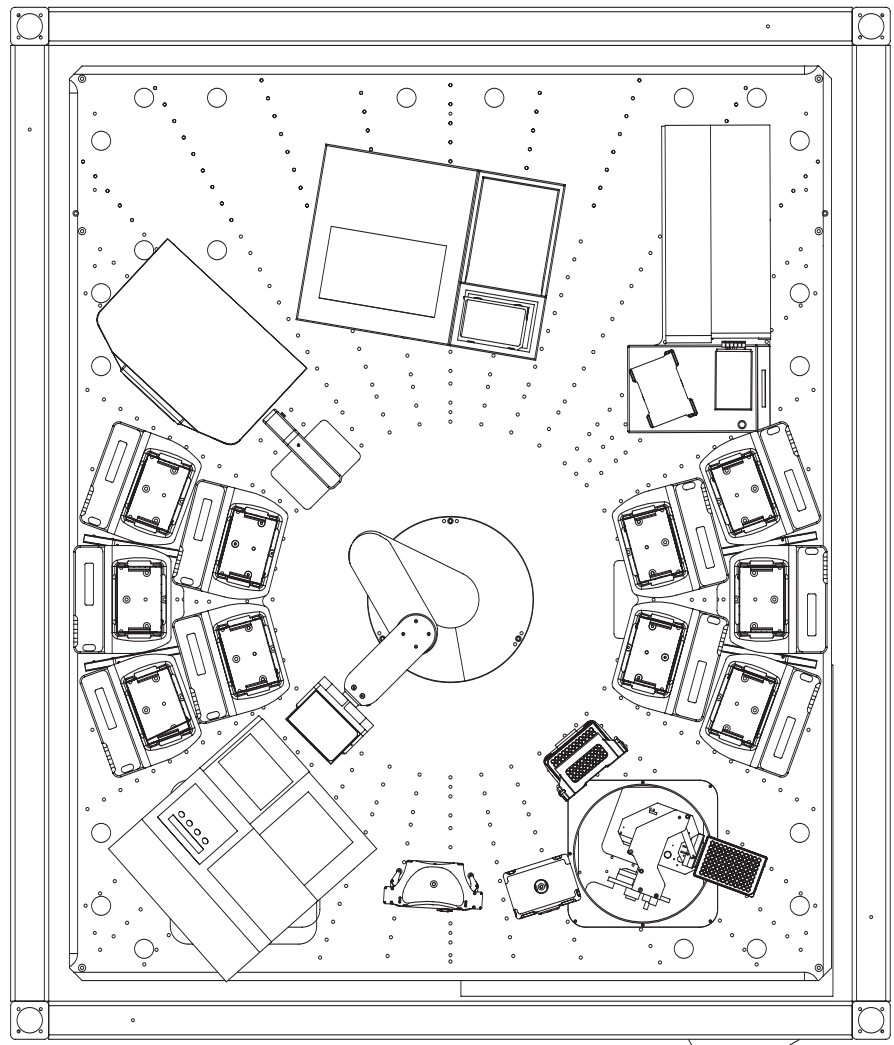


Agilent BioCel System Configuration

Cell-Based ELISA

Application Bulletin

The handling of cells in automated systems requires smooth and reliable operation and specialized equipment. Agilent Automation Solutions designed this BioCel to dispense cells into microplates and perform a series of alternating incubations and reagent additions. Intermediate plate batches, cell medium removal and absorbance detection is also part of the process. The BioCel is ideally suited to this complex assay format thanks to its dynamic scheduling software, flexible liquid handling options and the ability to integrate third-party instruments. In this particular application, the BioCel contains an Agilent Vertical Pipetting Station with tip-washing and auto-filling reservoir capabilities and 10 Agilent Labware Stacker units to accommodate large batches of plates. A plate washer (Bio-Tek Elx405), CO₂ incubator (LiCONiC StoreX) and a microplate absorbance reader (Thermo Multiskan Ascent) are integrated into the system. With the accompanying Agilent VWorks software and scheduler, this BioCel can handle approximately 500 plates per run in this 3-day experiment. Cells can be plated and stored at the rate of 1 min/plate.



Module List Component	Quantity	Function
Agilent robot	1	360° high speed robotic plate handler
Agilent Vertical Pipetting Station	1	Low volume sample transfer
Agilent Microplate Barcode Labeler	1	Plate identification and labeler
Bio-Tek ELx-405	1	Plate washer and bulk reagent dispenser
LiCONiC StoreX	1	240-plate storage
Agilent Labware Stacker	10	50-plate room-temperature storage
Thermo Multiskan Ascent	1	UV/visible absorbance reader
Lid Hotel*	1	Lid removal, storage and return

* Lid Hotel concept developed by Novartis Pharma AG, NIBR/DT/IAT, Basel, Switzerland

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.

Thermo and Multiskan Ascent are registered trademarks of Thermo. ELx405 is a trademark of Bio-Tek. StoreX is a trademark of Liconic Instruments.

© Agilent Technologies, Inc. 2009
Published in the U.S.A. February 25, 2009
5990-3624EN



Agilent Technologies