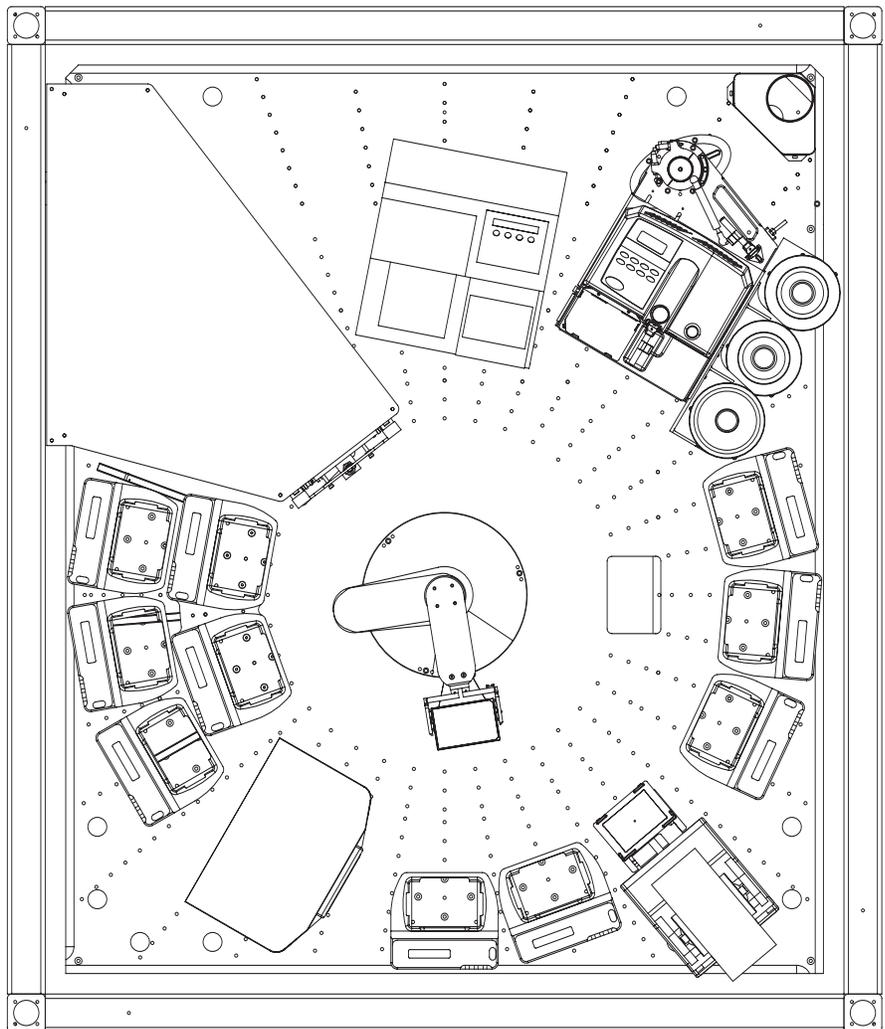


Agilent BioCel System Configuration

Flexible HTS System

Application Bulletin

The variety of different assay protocols for HTS is increasing constantly. Today's automated systems need to be flexible and cleverly designed to accommodate a large number of different protocols without major delays or costs involved. To prevent obsolescence, an automated system should be flexible enough to accommodate today's assays as well as new protocols and hardware. This Agilent BioCel includes a high quality selection of liquid handlers and readers, and an active environmental control for the entire system. A multimode reader (Zeiss Plate::Vision), bulk reagent dispensers (Genetix QFill2, Aliquot), plate washer (Bio-Tek ELx405 Select) and the Agilent Vertical Pipetting Station allow assays to be run in 96-, 384- or 1536-format. In combination with the multimode reader, a variety of detection modes are used: FI, FP, TRF and absorbance. The active environmental control allows assays to be run at constant temperature (ambient up to 37°C), avoiding edge effects often caused by transport into and out of incubators. The high precision, low volume pipettor Vertical Pipetting Station and the large plate storage capacity (Agilent Labware Stacker) allow the compound plate bank to be run alongside the assay, adding the compounds on-the-fly. No prior assay plate generation is necessary and due to the low volume capacity of the Vertical Pipetting Station (200 nL), DMSO concentrations in the assay are kept low.



Module List

Component	Quantity	Function
Agilent robot	1	360° high speed robotic plate handler
Agilent Vertical Pipetting Station	1	Low volume compound transfer
Genetix QFill2	2	Bulk reagent dispenser
Genetix aliquot	1	low volume bulk reagent dispenser
Agilent PlateLoc	1	Thermal plate sealer
Bio-Tek ELx405 Select	1	Plate washer
Vacuum Pad	1	Vacuum filtration station
Zeiss Multimode	1	Plate reader
Agilent Labware Stacker	10	50-plate storage
Environmental Control	1	Heated to 37°C, humidified to 75% r.h.

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.

Zeiss and Plate:Vision are registered trademarks of Carl Zeiss, Inc. ELx405 is a trademark of Bio-Tek.

© Agilent Technologies, Inc. 2009
Published in the U.S.A. February 25, 2009
5990-3619EN



Agilent Technologies