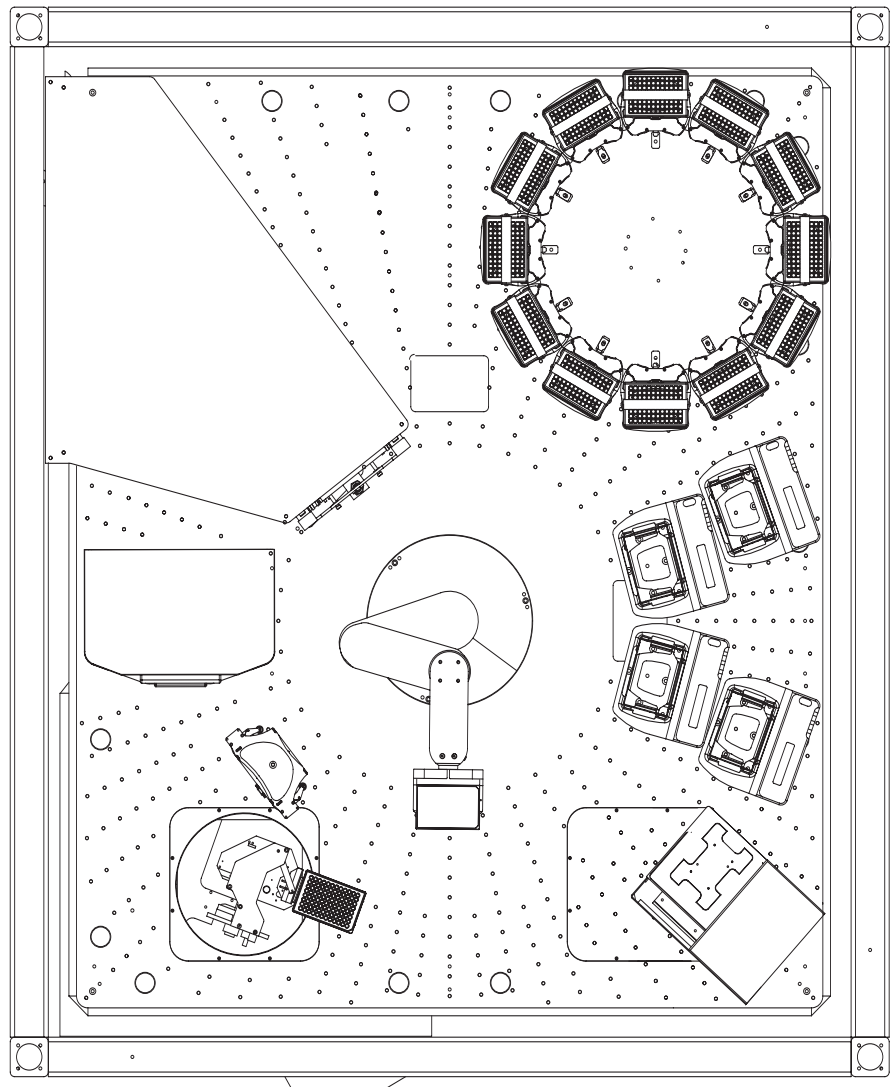


Agilent BioCel System Configuration Bioluminescent Kinase Assay Application Bulletin

There are a variety of kinase assay protocols utilized in laboratories today that address a variety of needs from high throughput screening to lead optimization. The flexibility of the Agilent BioCel System fits very nicely in this niche with its capability of changing and growing with the diverse needs of the assays and laboratories themselves.

Agilent Automation Solutions designed this BioCel System to perform a bioluminescent antibody-independent kinase assay. It utilizes an Agilent Vertical Pipetting Station to transfer compounds stored in 1536-well plates to 1536-well assay plates. The lids of the compound plates are removed and stored on the Lid Hotel as the plate is processed. During the temperature independent incubation steps, plates are moved into the PlateHub carousel and for incubation steps that require temperature control, they are moved into the Liconic incubator. Three Flexispense dispensers dispense kinase, substrate/ATP and stop/luciferase solutions, respectively. The emitted luminescence is quantified by the Zeiss Plate::Vision reader. The system can process a batch of 192 assay plates in 1536-well format in approximately 12 hours.



Module List

Component	Quantity	Function
Agilent robot	1	360° high speed robotic plate handler
Agilent Vertical Pipetting Station	1	Low volume liquid transfer
PlateHub	1	192-plate ambient storage
Lid Hotel*	1	Lid removal, storage and return
LiCONiC Incubator	1	240-plate controlled environment storage
Asys Hitech Flexispense	1	Low volume bulk reagent dispenser
Zeiss Plate::Vision	1	Multimode plate reader
Agilent Labware Stacker	4	50-plate capacity storage

* 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.

Zeiss and Plate::Vision are registered trademarks of Carl Zeiss, Inc.

© Agilent Technologies, Inc. 2009
Published in the U.S.A. February 25, 2009
5990-3620EN



Agilent Technologies