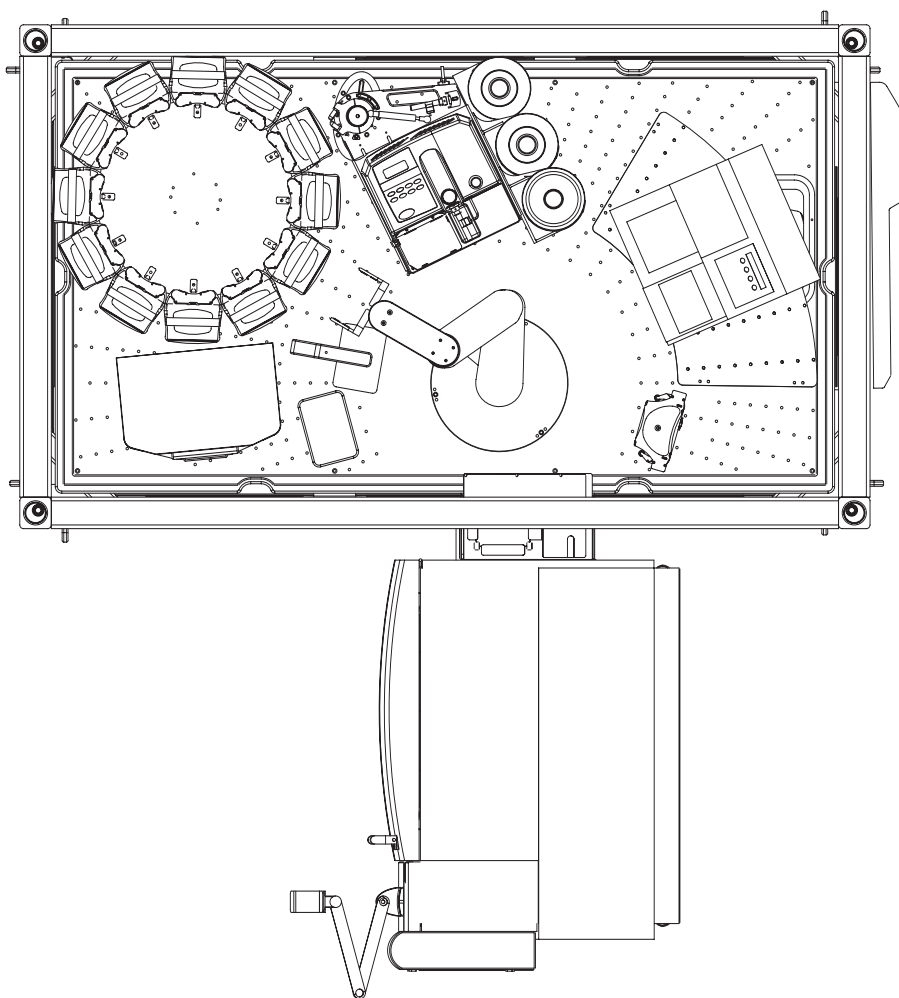


# Agilent BioCel System Configuration Automated FLIPR Assay in 384 Format Application Bulletin

The GPCR family is a very important target area for Drug Discovery and many of those targets can be assayed via intracellular calcium measurement. The FLIPR, and especially the new FLIPR<sup>Tetra</sup> from Molecular Devices is designed to perform this measurement in a reliable and effective manner. Agilent's BioCel System, with its speed and versatile scheduling software is a perfect match for the FLIPR<sup>Tetra</sup>, resulting in high quality output and very high throughput. The typical set-up of a calcium flux assay on the FLIPR can be run either in the classical way, including the addition of Fluo-4 and subsequent washing, or using the Molecular Devices Calcium-3 kit, omitting the washing steps. Both methods can easily be accommodated on this BioCel System. Addition of the loading buffer is handled by the bulk dispenser and the (optional) washing steps are executed with the Bio-Tek ELx405 plate washer. The Agilent Vertical Pipetting Station, with its high precision pipetting and low volume capacities, is ideally suited for compound additions, keeping DMSO concentrations in the assay low. The plate carousel and the Liconic incubator allow setting up assay batches of 192 plates to be completed in about 20 hours for a typical FLIPR antagonist screen.



## Module List

Component	Quantity	Function
Agilent robot	1	360° high speed robotic plate handler
Agilent Vertical Pipetting Station	1	Low volume compound transfer
Genetix aliQuot	1	Bulk reagent dispenser
Bio-Tek ELx405	1	Plate washer
FLIPR <sup>Tetra</sup>	1	Plate reader
PlateHub	1	192-plate capacity carousel
Liconic incubator	1	240-plate temp & CO <sub>2</sub> controlled incubation
Lid Hotel*	1	Lid removal, storage and replacement

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

[www.agilent.com/lifesciences/automation](http://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.

Molecular Devices, FLIPR, and FLIPR<sup>Tetra</sup> are registered trademarks of MDS Analytical Technologies. ELx405 is a trademark of Bio-Tek.

© Agilent Technologies, Inc. 2009  
Published in the U.S.A. February 25, 2009  
5990-3625EN



**Agilent Technologies**