

## Painting a picture of 20,000 genes in the mouse brain

Tecan's GenePaint™ systems for automated, high throughput *in situ* hybridization have been used to develop a groundbreaking, genome-wide atlas of gene expression in the adult mouse brain. The comprehensive digital atlas details the expression patterns of approximately 20,000 genes, providing an invaluable source of information for scientists studying brain organization, gene functions and behavior. Watch out for a special feature on the project in the next issue of the Tecan Journal!



## 7 detection modes

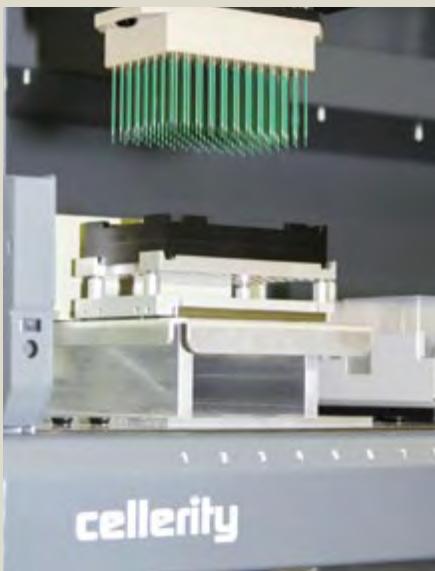
Tecan's highly sensitive and extremely fast Safire<sup>2</sup>™ and Infinite™ F500 microplate readers offer seven different detection modes, saving valuable time for a wide range of high throughput applications in the biotech, pharmaceutical, research and cosmetics industries. These two modular and upgradeable systems provide very reliable performance with top quality results and are easily integrated into Tecan's Freedom EVO® series of liquid handling workstations.



Relax  
your reader is a Tecan

[www.tecan.com/relax](http://www.tecan.com/relax)

## Cellerity™ is now available with the MultiChannel Arm™



The Cellerity™, Tecan's fully automated cell culture system, is now available with the MultiChannel Arm™ (MCA) module, increasing the speed and throughput of plating cells into microplates for assays. The MCA module has a 96-channel pipetting head and can be operated with washable and disposable tips. The new option is compatible with lidded and insert plates, and all standard Cellerity options, delivering a greater level of flexibility to your cell culture platform.

