

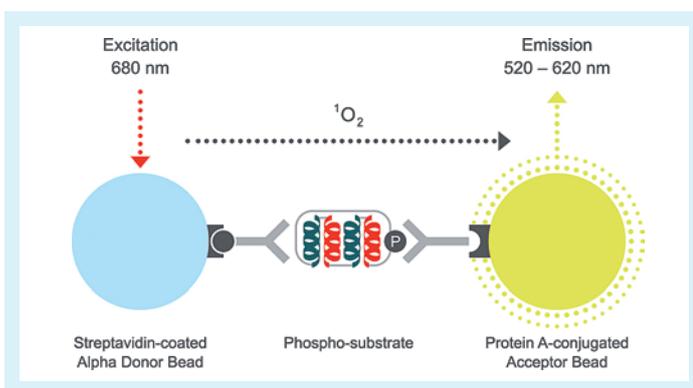
# AlphaScreen® option launched for Infinite F200 PRO multimode reader

The newly developed AlphaScreen module for the popular Infinite F200 PRO multimode reader has been designed to provide customers with a cost-effective solution for AlphaScreen and AlphaLISA® assays. Offering excellent sensitivity for screening and research, the new option further extends the flexibility of the Infinite F200 PRO for low and medium throughput applications.

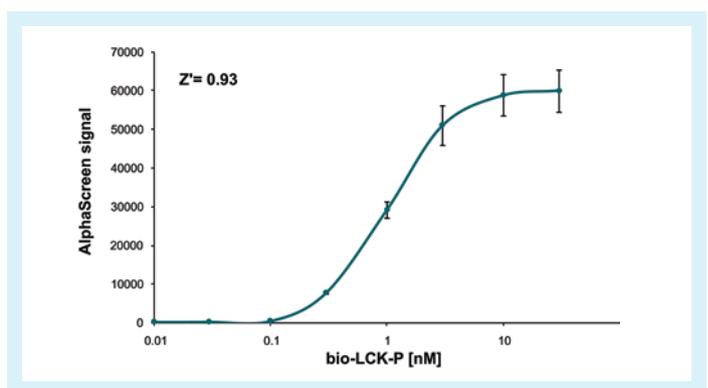
The AlphaScreen module combines the Infinite F200 PRO's established fluorescence top optics with a bespoke filter set designed to use the instrument's standard flashlamp. This set-up ensures optimal assay performance without the need for a specialized reader, complementing the

system's absorbance, fluorescence intensity, luminescence, fluorescence polarization and time-resolved fluorescence resonance energy transfer measurement modes.

To find out more on Tecan's AlphaScreen solutions, visit [www.tecan.com/alphascreen](http://www.tecan.com/alphascreen)



The bead-based AlphaScreen assay system



A typical AlphaScreen signal curve using an Infinite F200 PRO with PerkinElmer's P-Tyr-100 phosphotyrosine kinase detection kit

## First Air LiHa modules shipped to customers

Tecan is pleased to announce that the first Air LiHa air displacement pipetting arms have left Tecan's production facility in Männedorf. Designed to offer an unrivalled level of flexibility for Freedom EVO liquid handling workstations, this innovative module is Tecan's most compact single channel pipetting option, and offered a number of fresh challenges for the manufacturing team. Kaspar Städeli, Assembly Team Leader, commented: "The Air LiHa is a very compact unit. Like a good Swiss watch, many small parts need to work in perfect unison, and we have invested many hours in training and in discussion with our trusted suppliers to guarantee the highest quality standards. We have created a whole new generation of manufacturing tools and testing equipment for this project, furthering our precision engineering capabilities and ensuring that the Air LiHa offers low maintenance, high performance operation."



The Air LiHa Assembly Team (back row, left to right: Sven Schmidt, Claudio Monti, Firat Sert, Kaspar Städeli, Thorsten Jörg, Jörg Egli) together with Project Leader Urs Schumacher (front left) and Product Manager Ralf Masantschek (front right)

Fredy Hottinger, Head of Assembly in Männedorf, added: "The transfer of any new product from development to production is always a critical step in terms of product quality. We have worked more closely with R&D than ever before on the Air LiHa – it has been a real

team effort – and we are proud of the quality and performance of this exciting new offering, and the benefits it will bring to our customers."

To find out more on Tecan's Air LiHa, visit [www.tecan.com/airliha](http://www.tecan.com/airliha)