

Tecan Microplates.

ASSURED PERFORMANCE FOR ABSORBANCE, FLUORESCENCE AND LUMINESCENCE MEASUREMENTS - PLUS CELL IMAGING - IN ANSI/SLAS FORMAT.



Tecan Microplates.

Tecan Microplates are the latest addition to the Tecan consumables portfolio. We offer a selection of polystyrene, medium-binding ANSI/SLAS-format microplates compatible with Tecan instruments.

- Optimal plate height and manufacturing tolerances allow reader optics to be as close to the plate as possible, avoiding well-to-well signal crosstalk.
- Developed and tested to work in combination with the imaging algorithms of our multimode readers, offering assured performance.
- Microplate well diameter is optimal to work with the Spark® system's bright field imaging module - critical for confluence assessments.

ENSURE RELIABILITY OF YOUR DATA BY USING TECAN MICROPLATES - COMPATIBLE WITH THE SPARK MULTIMODE READER.

Obtaining as much data as possible is essential to ensure comprehensive results for automated confluence measurements. Microplate well diameter plays an important role in this, and so Tecan Microplates are optimized to maximize the number of data points and images that can be captured with the Spark platform (see Figure 1).

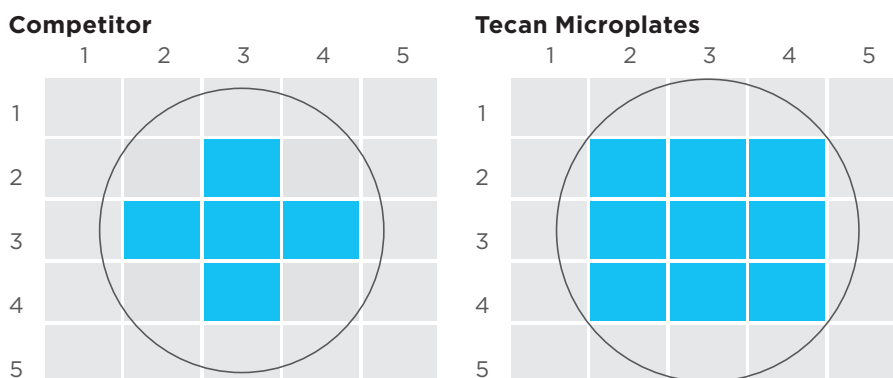


Figure 1: Plate well diameter is critical for accurate analysis.

Biochemical Assay Plates.

Black and white biochemical assay plates are designed for fluorescence and luminescence measurement assays.

The low auto-phosphorescence of our white biochemical plates helps to reduce pre-measurement incubation times, while our unique combination of black microplates plates and black lids allows uninterrupted measurement of light-sensitive assays using Spark's unique Lid Lifter™ system.

WHITE BIOCHEMICAL ASSAY PLATES

Most white microplates load up with photons when kept under normal laboratory light. This can critically influence the background of luminescence assays, especially when measured in highly sensitive luminescence reader, such as Spark.

Designed for luminescence, TR-FRET and Alpha technology assays, Tecan white biochemical assay plates enable immediate measurements while still offering good results (see Figure 2).

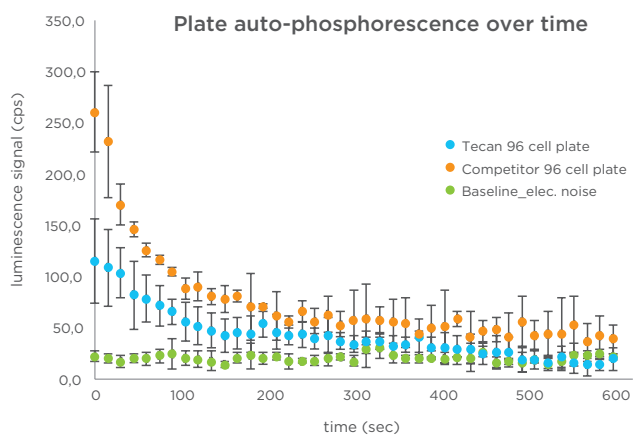

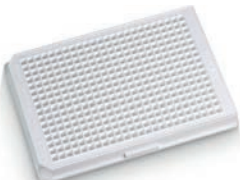


Figure 2: Low auto-phosphorescence of white Tecan Microplates compared to a competitor plate.



Product selection

	Product	Product
	96-well, white, flat bottom 10 pcs/bag, 40 pcs/box Material number: 30122300	
		384-well, white, flat bottom 10 pcs/bag, 40 pcs/box Material number: 30122301

BLACK BIOCHEMICAL ASSAY PLATES

Intended for use in fluorescence assays, including fluorescence polarization (FP), time-resolved fluorescence (TRF) and TR-FRET applications.

Product selection

	Product		Product
	96-well, black, flat bottom 10 pcs/bag, 40 pcs/box Material number: 30122298		384-well, black, flat bottom 10 pcs/bag, 40 pcs/box Material number: 30122299

BIOCHEMICAL ASSAY PLATE LIDS

Tecan biochemical assay plates are available with a choice of two low-profile plate lids – black and clear – helping to ensure that assays with long incubation times will not be affected by evaporation.

We offer a convenient solution for fluorescence assays that require long incubations in the dark. Simply use a black biochemical assay plate and lid in combination with Spark's Lid Lifter, and incubate for as long as required. Using a magnetic disc attached to the plate lid, the Lid Lifting can automatically remove the plate lid to perform measurements at specified time intervals.

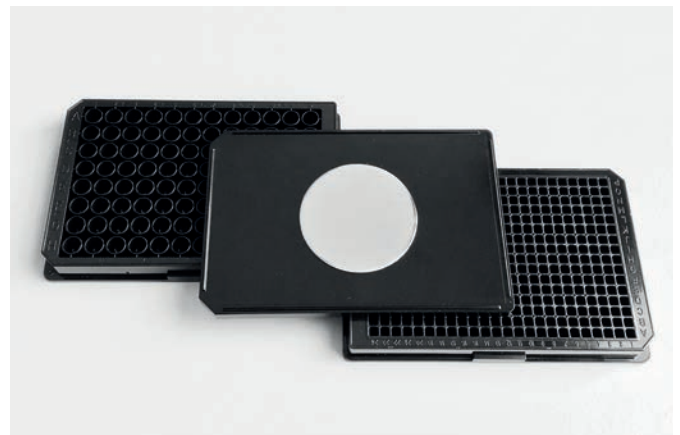





Figure 3: Black Tecan Microplates and the black Tecan lid with the lid lifting disc attached.

Product selection

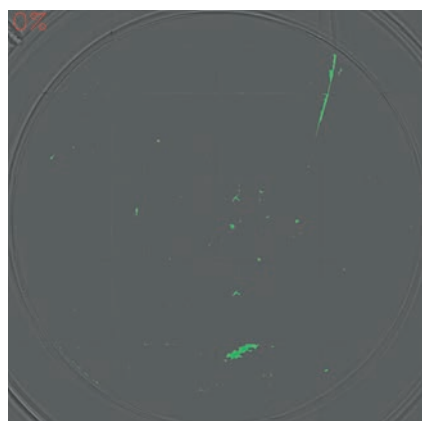
	Product		Product
	Biochemical assay plate lids, clear 20 pcs/bag, 80 pcs/box Material number: 30122308		Lid Lifter discs 50 pcs/box Material number: 30090494
	Biochemical assay plate lids, black 20 pcs/bag, 80 pcs/box Material number: 30122309		

Cell Culture Plates.

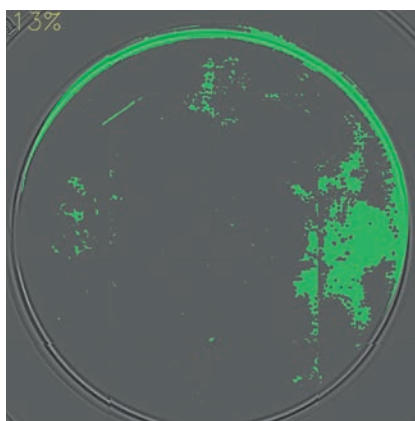
A collection of tissue culture-treated (TC-treated), sterile cell culture plates with individual lids for confluence and bottom-reading fluorescence measurements in cell-based applications.

CLEAR CELL CULTURE PLATES

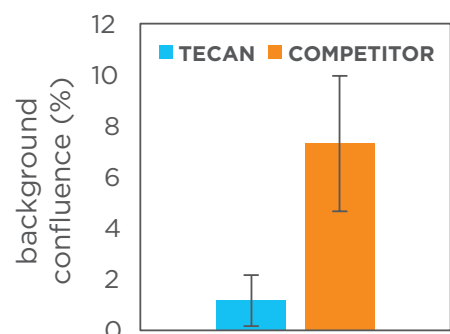
It is critical to ensure the optical quality of microplate well bottoms for contrast-based confluence measurements, as optical artifacts or particles that have a higher contrast than the background could be incorrectly identified as cells. Our cell culture plates offer assured optical performance, with minimum background signal.



Tecan Microplate







Competitor C Microplate



	Tecan	Competitor C
Average	1.2	7.3
SD	0.9	2.6


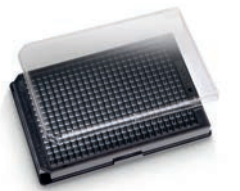
Figure 4: Difference in background confluence (artifacts) between a Tecan cell culture plate and a competitor plate using the Spark reader.

Product selection

	Product		Product
	24-well, clear with lid, flat bottom, TC-treated, sterile Individually packed, 100 pcs/box Material number: 30122302		96-well, clear with lid, flat bottom, TC-treated, sterile Individually packed, 100 pcs/box Material number: 30122304
	48-well, clear with lid, flat bottom, TC-treated, sterile Individually packed, 100 pcs/box Material number: 30122303		384-well, clear with lid, flat bottom, TC-treated, sterile 8 pcs/bag, 32 pcs/box Material number: 30122305

BLACK CELL CULTURE PLATES WITH CLEAR FILM BOTTOM

Product selection

	Product		Product
	96-well, black with lid, clear flat bottom, TC-treated, sterile 8 pcs/bag, 32 pcs/box Material number: 30122306		384-well, black with lid, clear flat bottom, TC-treated, sterile 8 pcs/bag, 32 pcs/box Material number: 30122307



We are offering a 20% discount on your first order!

Tecan Multimode Readers.

Infinite® F50

The Infinite F50 and Infinite F50 Robotic are Tecan's family of reliable, validated and compact absorbance microplate readers for ELISA. They provide accurate, fast photometric measurements using innovative eight-channel optics and LED technology.



Sunrise™

Tecan's Sunrise absorbance microplate reader offers all the functionality needed for numerous photometric applications, including advanced 12-channel optics for fast, multichannel absorbance reading of ELISAs, a temperature control function for enzyme kinetic assays, and a tunable wavelength function for wavelength scanning.



Infinite® M Plex

The Infinite 200 PRO is an easy-to-use multimode plate reader family that offers affordable, high performance detection solutions, referenced in more than 1800 peer reviewed publications.



Spark®

The Spark multimode microplate reader is your research partner – built for today and ready for tomorrow. It can effortlessly handle your routine or complex assays, with a modular and fully upgradeable design. Its unique technology and software features ensure that there is a Spark configuration to increase the productivity of your lab, both today and in the future.



Contact your local Tecan representative to order Tecan Microplates or request a quote.
For more details, visit the Tecan Consumables webpage: www.tecan.com/consumables

.....
Australia +61 3 9647 4100 **Austria** +43 62 46 89 330 **Belgium** +32 15 42 13 19 **China** +86 21 220 63 206 **France** +33 4 72 76 04 80 **Germany** +49 79 51 94 170
Italy +39 02 92 44 790 **Japan** +81 44 556 73 11 **Netherlands** +31 18 34 48 17 4 **Nordic** +46 8 750 39 40 **Singapore** +65 644 41 886 **Spain** +34 93 595 25 31
Switzerland +41 44 922 89 22 **UK** +44 118 9300 300 **USA** +1 919 361 5200 **Other countries** +41 44 922 81 11
.....

Tecan Group Ltd. makes every effort to include accurate and up-to-date information within this publication, however, it is possible that omissions or errors might have occurred. Tecan Group Ltd. cannot, therefore, make any representations or warranties, expressed or implied, as to the accuracy or completeness of the information provided in this publication. Changes in this publication can be made at any time without notice. All mentioned trademarks are protected by law. In general, the trademarks and designs referenced herein are trademarks, or registered trademarks, of Tecan Group Ltd., Männedorf, Switzerland. A complete list may be found at <http://www.tecan.com/trademarks>. Product names and company names that are not contained in the list but are noted herein may be the trademarks of their respective owners. For technical details and detailed procedures of the specifications provided in this document please contact your Tecan representative.

Tecan is in major countries a registered trademark of Tecan Group Ltd., Männedorf, Switzerland.
© 2017 Tecan Trading AG, Switzerland, all rights reserved.

www.tecan.com



400217 V2.0 06-2017