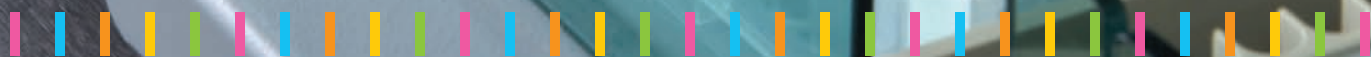




# Tecan AC

# Extraction Plate™

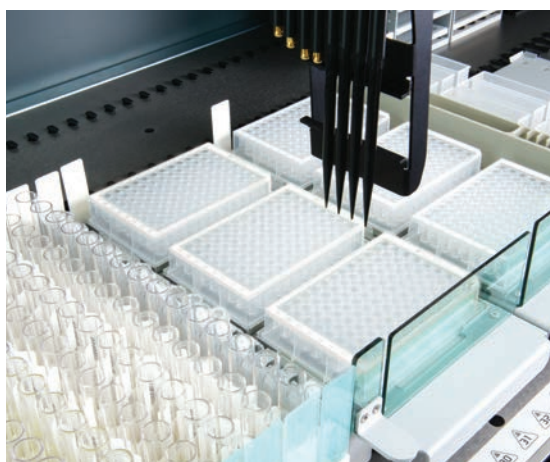
SAMPLE PREPARATION FOR LC-MS MADE EASY



## AC Extraction Plate for higher throughput, faster sample preparation

The AC Extraction Plate is a convenient, automation-friendly solution for your LC-MS sample preparation challenges. It reduces the process to a simple 'pipette and shake' routine, eliminating the need for filtration, centrifugation and solvent evaporation steps.

### Sample preparation in 6 easy steps



Thanks to its compact standard ANSI/SLAS format, the AC Extraction Plate does not impact your workbench design and provides you with a flexible choice of carrier positions.



Method development and protocol optimization for a new analyte can be easily done manually before transferring the process to the automated liquid handler.

### KEY BENEFITS

#### Flexible method development

The AC Extraction Plate is suitable for both automated and manual processes. Transferring your manually optimized protocols to an automated liquid handling platform has never been easier.

#### Ideal for high throughput

The standard ANSI/SLAS format of the AC Extraction Plate allows easy automation on a liquid handling platform such as the Freedom EVO®.

#### Convenient and fast

After unpacking, the AC Extraction Plate can be used immediately, without the need for conditioning or other pre-treatments.

#### A wide range of applications

The AC Extraction Plate offers an open development platform that can be applied to sample preparation for a broad range of small molecule analytes from biological fluids such as vitamin D, steroids, immunosuppressants and other small molecules.

#### Faster and easier workflow

The sample preparation routine consists of simple pipette and shake steps, all performed in one microplate. This eliminates the need for sample transfers, centrifuges or vacuum systems.

#### Highly reproducible results

Sample preparation is easily standardized and automated with the AC Extraction Plate to eliminate random errors and increase consistency of results.

## How does the AC Extraction Plate based on TICE™ technology work?

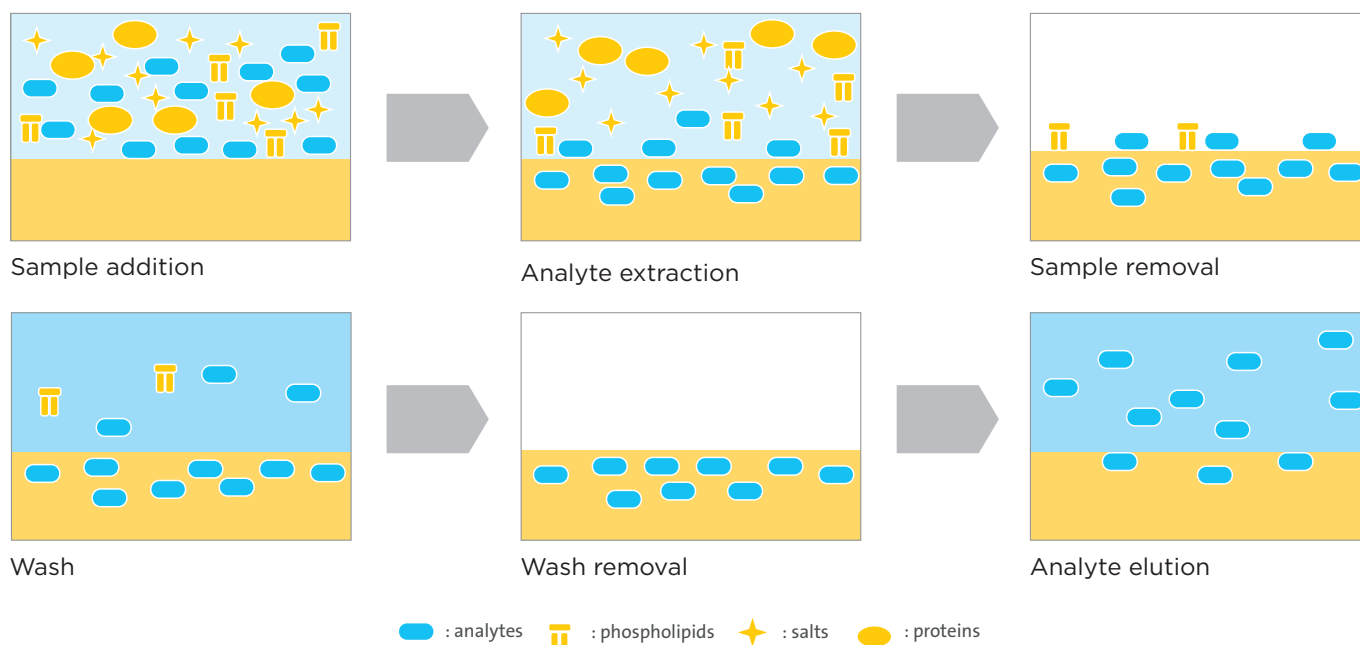
The AC Extraction Plate is a convenient, automation-friendly solution for your LC-MS sample preparation challenges. It reduces the process to a simple 'pipette and shake' routine, eliminating the need for filtration, centrifugation and solvent evaporation steps.

### Sample preparation process

Sample preparation is a straightforward three-step process. First, the analyte is **extracted** from a biological fluid, such as serum, plasma, urine or whole blood.

This is followed by a **wash** step to ensure removal of matrix materials – such as proteins, phospholipids and salts.

And, finally, **elution** from the absorptive coating by a specifically selected solvent mixture. The resulting eluate is of sufficiently high purity that it can be injected directly into an LC-MS system for analysis.



Tecan AC Extraction Plate.



Graphical representation of the Tecan AC Extraction Plate (left) and a cross section of a single well (right), showing the presence of the TICE coating to approximately half the well's height.

## SPECIFICATIONS

<b>Chemical resistance</b>	Known compatible solvents: methanol, acetonitrile, ethanol, isopropanol, 2% acetic acid 2% formic acid, 2% ammonium hydroxide, ammonium acetate Known incompatible solvents: aromatic hydrocarbons, esters, ketones and concentrated acids
<b>Operational conditions</b>	Standard laboratory range 15° to 32°C (59° to 90°F). The relative operating humidity must be 30 to 80 % (non-condensing).
Storage:	15° to 32°C (59° to 90°F). Must be kept dry and protected against direct sunlight.
Tested pH range:	Aqueous solutions with a pH range of 2 to 10
Max. liquid volume:	500 µl
Expiration date:	The AC Extraction Plate can be used for up to 12 months from the date of manufacture when properly stored. The expiration date is indicated on the primary packaging label.
Number of tests:	96 tests
Dimensions:	127.8 x 85.5 x 27.1 mm (WxDxH)
<b>Order information</b>	
Material number:	30072211
Package contents:	AC Extraction Plate, 96-wells with TICE coating, 1 plate per unit, Certificate of Analysis, instructions for use.



The AC Extraction Plate is sealed in an aluminum protective foil and packaged in a robust box together with the Certificate of Analysis and instructions for use.

## Tecan – Who we are

Tecan is a leading global provider of laboratory instruments and solutions in biopharmaceuticals, forensics, clinical diagnostics, academic centers and life science industries, and specializes in the development and production of automation solutions, detection instruments such as microplate readers, microarray-related products and washers.

Founded in Switzerland in 1980, Tecan has manufacturing, research and development sites in both North America and Europe, and maintains a sales and service network in 52 countries. To date, Tecan has distributed approximately 20,000 microplate readers worldwide and is committed to continuous technological improvements and compliance to global quality standards.

.....  
**Australia** +61 3 9647 4100 **Austria** +43 62 46 89 33 **Belgium** +32 15 42 13 19 **China** +86 21 220 63 206 **Denmark** +45 70 23 44 50 **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 **Singapore** +65 644 41 886 **Spain** +34 93 490 01 74  
**Sweden** +46 8 750 39 40 **Switzerland** +41 44 922 89 22 **UK** +44 118 9300 300 **USA** +1 919 361 5200 **Other countries** +43 62 46 89 33  
.....

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. For technical details and detailed procedures of the specifications provided in this document please contact your Tecan representative. This brochure may contain reference to applications and products which are not available in all markets. Please check with your local sales representative. 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 [www.tecan.com/trademarks](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.

© 2017, Tecan Trading AG, Switzerland, all rights reserved. For disclaimer and trademarks please visit [www.tecan.com](http://www.tecan.com)

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

The Tecan logo consists of a red circle with a white dot inside, followed by the word "TECAN" in a bold, black, sans-serif font.

