

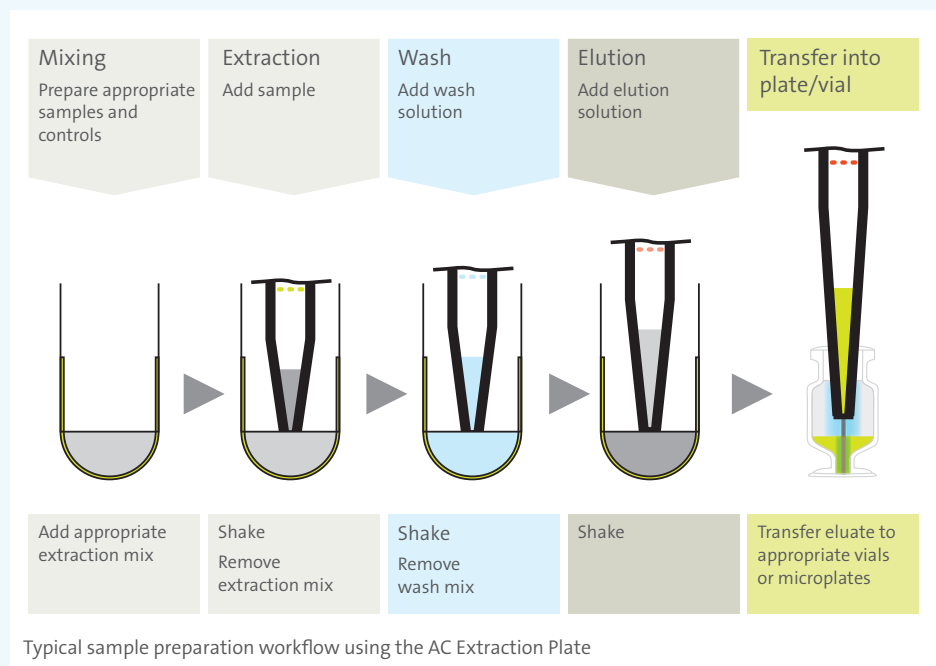
AC Extraction Plate™ – a unique sample preparation tool for LC-MS



The AC Extraction Plate offers straightforward sample preparation for LC-MS

The recently launched AC Extraction Plate – based on Tecan's proprietary TICE™ (Tecan Immobilized Coating Extraction) technology – is a convenient, automation-friendly solution designed to streamline sample preparation for LC-MS analysis of small molecules. The plate reduces the sample preparation process to a simple 'pipette and shake' routine, eliminating the need for laborious filtration, centrifugation and solvent evaporation steps.

The AC Extraction Plate is a 96-position, deep-well microplate, with the inner surface of each well coated with a highly controlled layer of TICE material. This coating efficiently extracts low molecular weight analytes – such as vitamin D and testosterone – from aqueous solutions and eliminates the need for protein precipitations and liquid-liquid separations. By avoiding these time-consuming and difficult to automate processes, the AC Extraction Plate provides a convenient and easy to automate solution for the preparation of biological samples and other complex fluids for LC-MS analysis of small molecules.



To find out more about Tecan's AC Extraction Plate, visit www.tecan.com/acplate