

Extraction of hydrophobic bases from complex liquid samples with SOLA WCX SPE

Author

Thermo Fisher Scientific

Keywords

Solid Phase Extraction (SPE), small molecule purification, extraction of hydrophobic bases

Introduction

Thermo Scientific™ SOLA™ is a solid-phase extraction (SPE) cartridge featuring mixed-mode polymeric sorbent and a fritless design for small sample sizes (typically 25–400 μL). It is available in SOLA 10 mg and Thermo Scientific™ SOLA μ ™ 2 mg formats. The fritless design reduces hold-up volume and improves consistency of extraction. The SOLA WCX has reversed-phase (RP) and weak cation exchange (WCX) functions. The typical use is for the extraction of hydrophobic bases from complex liquid samples.

Important notes

- Maximum loading capacity is ~10% of sorbent weight
- Sample should be processed through the cartridge at about 1 mL/min; too high a flow can lead to inconsistent results
- The volumes given are typical, and should be optimized for the analyte and matrix of interest

Materials required

- Methanol, LCMS grade
- 2% formic acid in methanol, LCMS grade
- 5% ammonium hydroxide in water, LCMS grade
- 10–30% acetonitrile in water (optional), LCMS grade
- SPE vacuum manifold, vacuum regulator, vacuum pump
- 96-well collection plate, appropriate to final extract volume, 25–200 μL per sample
- Waste-collection tray or plate, ~1800 μL per sample
- Pipettes and tips

Protocol

1. Prepare the sample for extraction
 - Dilute viscous samples (e.g., plasma) 1:1 with water
 - When the analytes bind to matrix proteins, 1:1 dilution with 10–30% aqueous acetonitrile can improve recovery
 - Adjust to a $\text{pH} \leq \text{pKa} - 2$ and $\text{pH} \geq 7$ as necessary to ionize the analytes and SPE
 - Add internal standard if desired
2. Prepare the SOLA SPE for sample loading
 - Wash with $2 \times 100 \mu\text{L}$ of methanol (optional)
 - Wash with $2 \times 100 \mu\text{L}$ of 5% ammonium hydroxide in water. Do not let cartridge dry before loading sample.
3. Load the sample onto the SOLA SPE at a flowrate of about 1 mL/min
 - Up to 800 μL of prepared sample
4. Wash away interferences
 - Wash with $2 \times 100 \mu\text{L}$ of 5% ammonium hydroxide in water. This removes salts, acids, proteins, carbohydrates.
 - Wash with $2 \times 100 \mu\text{L}$ of methanol. This removes hydrophobic, neutral and anionic matrix components. Alkaline methanol can improve recovery of cationic analytes for some cases. Let cartridge dry a few minutes before elution.
5. Collect analyte fraction in the sample well plate
 - Elute with $\geq 2 \times 12.5 \mu\text{L}$ (SOLA μ) or $\geq 2 \times 50 \mu\text{L}$ (SOLA) of 2% formic acid in methanol. Elute each aliquot initially by gravity then apply vacuum/pressure to ensure all solvent is eluted from the cartridge.
6. Post-extraction
 - If necessary, evaporate and re-constitute in a compatible solvent
 - For RP-LC analysis, dilute to $\leq 50\%$ organic solvent

Related products

Description	Part Number
SOLA WCX, 10 mg/1 mL Cartridges	60109-004
SOLA μ WCX, 10 mg/1 mL 96 Well Plates	60209-004
SOLA WCX, 10 mg/2 mL 96 Well Plates	60309-004
Thermo Scientific™ Hypersep™ Universal SPE Vacuum Manifold, for 96-well plate or 24/48 cartridges	60104-230
Thermo Scientific™ Hypersep™ Vacuum Pump, European version	60104-241
Thermo Scientific™ Hypersep™ Vacuum Pump, North American version	60104-243

Current versions of product instructions are available at separatedbyexperience.com/chromexpert

Learn more about SOLA and SOLA μ Solid Phase Extraction at thermofisher.com/solaspe

ThermoFisher
SCIENTIFIC