thermoscientific

APPLICATION BRIEF 21933

Extraction of hydrophobic acids from complex liquid samples with SOLA WAX SPE

Author

Thermo Fisher Scientific

Keywords

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

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[™] SOLAµ[™] 2 mg formats. The fritless design reduces hold-up volume and improves consistency of extraction. The SOLA WAX has reversed-phase (RP) and weak anion exchange (WAX) functions. The typical use is for the extraction of hydrophobic acids 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 water, LCMS grade
- 5% ammonium hydroxide in methanol, 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



thermo scientific

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 pH ≥ pKa + 2 and pH < 9 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 μL of methanol (optional)
 - Wash with 2 \times 100 μL of 2% formic acid. Do not let cartridge dry before loading sample
- 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 μL of 2% formic acid in water. This removes salts, bases, proteins, carbohydrates.
 - Wash with 2 × 100 μL of methanol. This removes hydrophobic, neutral and cationic matrix components. Acidified methanol can improve recovery of anionic analytes for some cases. Let cartridge dry a few minutes before elution.
- 5. Collect analyte fraction in the sample well plate
 - Elute with ≥2 × 12.5 µL (SOLAµ) or ≥2 × 50 µL (SOLA) of 5% ammonium hydroxide 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 ≤50% organic solvent

Related products

Description	Part Number
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 SOLAµ Solid Phase Extraction at **thermofisher.com/solaspe**



© 2019 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries. This information is presented as an example of the capabilities of Thermo Fisher Scientific products. It is not intended to encourage use of these products in any manners that might infringe the intellectual property rights of others. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representatives for details. **AB21933-EN 0519M**