thermoscientific

SOLA and SOLAµ SPE method development

Consistent excellence for bioanalysis



Thermo Scientific™ SOLA™ products are designed for bioanalytical and clinical research analysts who are tasked with providing high-quality analytical results from complex biological samples in a high-throughput environment, while complying with strict legislation. These demands are compounded by the continued push to higher efficacy drugs and long acting formulations which continue to drive sensitive requirements to lower levels to enable accurate quantification.



In order to meet these demands, bioanalytical methods must provide:

SOLA SCY

- Robustness low analytical failure rates
- Ability to process low sample volumes
- High-sensitivity
- High-reproducibility
- Ease-of-use
- High-throughput processing
- Efficient and fast processes

Acidic co	Acidic compounds		Basic compounds		compounds Basic compounds	
Strong acids	Weak acids	Neutral compounds	Weak bases	Strong bases		
$(pK_a < 2)$	(pK _a 2-4)		(pK _a 8-10)	(pK _a > 10)		

The SOLA solid phase extraction (SPE) range meets these demands due to the unique and innovative frit-less SPE technology which eliminates the issues with traditional loose-packed SPE formats (see figure to the right). Combining the support material and active media components into a solid, uniform sorbent bed provides stable and controllable flow through characteristics and has an added advantage when dealing with viscous biological samples, as it prevents blocking and enables high throughput processing.

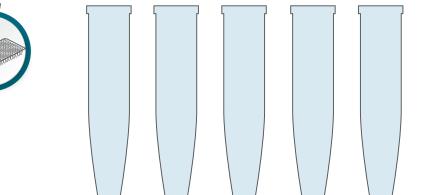
The manufacturing process also allows for high levels of reproducibility, not only from cartridge to cartridge or well to well but also batch to batch. The Thermo Scientific™ SOLAµ™ SPE range has the added benefit of being able to provide:

SOLA SAY

• Up to a 20 fold increase in sensitivity

SOLA WAY

- Ability to process samples restricted in volume
- Increased workflow efficiency and sample integrity



lucible positioning me after time	No voiding	No channeling	Reproducible packing time after time

SOLA WCY

SOLA WAX	SULA SAX	SOLA HRP	SOLA SCX	SOLA WCX				
Mixed-mode weak anion exchange	Mixed-mode strong anion exchange	Reversed-phase	Mixed-mode strong cation exchange	Mixed-mode weak cation exchange				
	Generic method protocol for cartridge and 96 well plate formats							
Condition 500 µL methanol	Condition 500 μL methanol	Condition 500 µL methanol	Condition 500 μL methanol	Condition 500 µL methanol				
Equilibrate 500 μL water with 1% formic acid	Equilibrate 500 μL water with 1% ammonium hydroxide	Equilibrate 500 μL water	Equilibrate 500 μL water with 1% formic acid	Equilibrate 500 µL water with 1% ammonium hydroxide				
Load 50-500 μL of sample at 1 mL/min	Load 50-500 μL of sample at 1 mL/min with 1% ammonium hydroxide	Load 50-500 μL of sample at 1 mL/min	Load 50-500 μL of sample at 1 mL/min with 1% formic acid	Load 50-500 µL of sample at 1 mL/min with 1% ammonium hydroxide				
Wash 1 500 μL water	Wash 1 500 μL water with 1% ammonium hydroxide	Wash 1 500 µL 5% methanol in water	Wash 1 500 µL water with 1% formic acid	Wash 1 500 μL water with 1% ammonium hydroxide				
Wash 2 500 μL methanol	Wash 2 500 μL methanol with 1% ammonium hydroxide		Wash 2 500 μL methanol with 1% formic acid	Wash 2 500 μL methanol with 1% ammonium hydroxide				
Elute 200-500 µL methanol with 1% ammonium hydroxide	Elute 200-500 μL methanol with 1% formic acid	Elute 200-500 μL methanol	Elute 200-500 µL methanol with 1% ammonium hydroxide	Elute 200-500 μL methanol with 1% formic acid				

SOLA HRD

SOLAµ SPE 96 Well Plates

Description	Bed weight	Well volume	Cat. no.	Quantity
SOLAµ HRP	2 mg	1 mL	60209-001	1 each
SOLAµ SCX	2 mg	1 mL	60209-002	1 each
SOLAµ SAX	2 mg	1 mL	60209-003	1 each
SOLAµ WCX	2 mg	1 mL	60209-004	1 each
SOLAµ WAX	2 mg	1 mL	60209-005	1 each

SOLA SPE 96 Well Plates

Description	Bed weight	Well volume	Cat. no.	Quantity
SOLA HRP	10 mg	2 mL	60309-001	1 each
SOLA SCX	10 mg	2 mL	60309-002	1 each
SOLA SAX	10 mg	2 mL	60309-003	1 each
SOLA WCX	10 mg	2 mL	60309-004	1 each
SOLA WAX	10 mg	2 mL	60309-005	1 each

SOLA SPE Cartridges

301/10: 1 Gai ti lagos					
Description	Bed weight	Well volume	Cat. no.	Quantity	
SOLAµ HRP	10 mg	1 mL	60109-001	100 pack	
SOLAµ SCX	10 mg	1 mL	60109-002	100 pack	
SOLAµ SAX	10 mg	1 mL	60109-003	100 pack	
SOLAµ WCX	10 mg	1 mL	60109-004	100 pack	
SOLAµ WAX	10 mg	1 mL	60109-005	100 pack	

