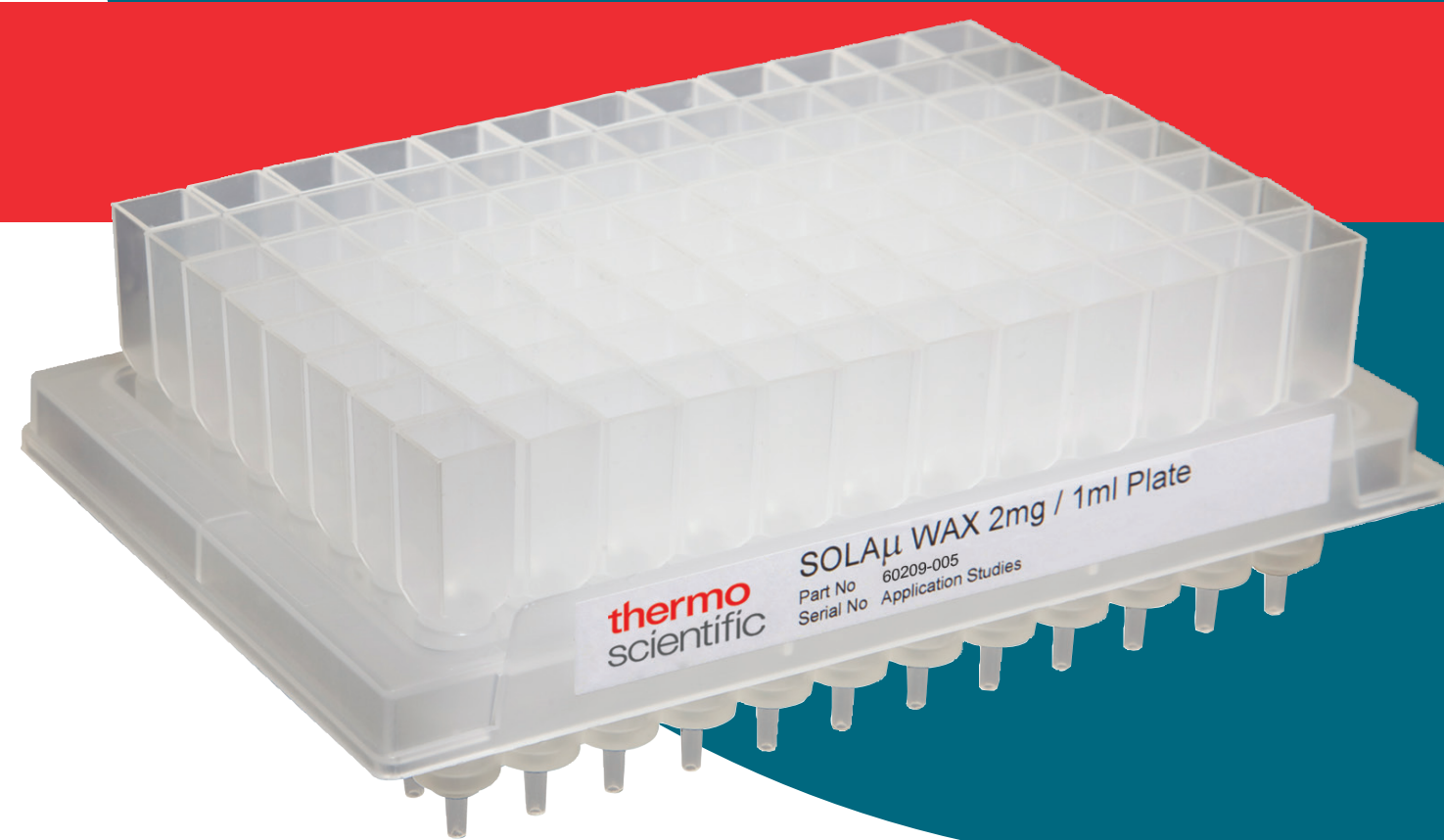


# SOLA and SOLA $\mu$ SPE method development

Consistent excellence for bioanalysis



## SOLA $\mu$ well plate

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:

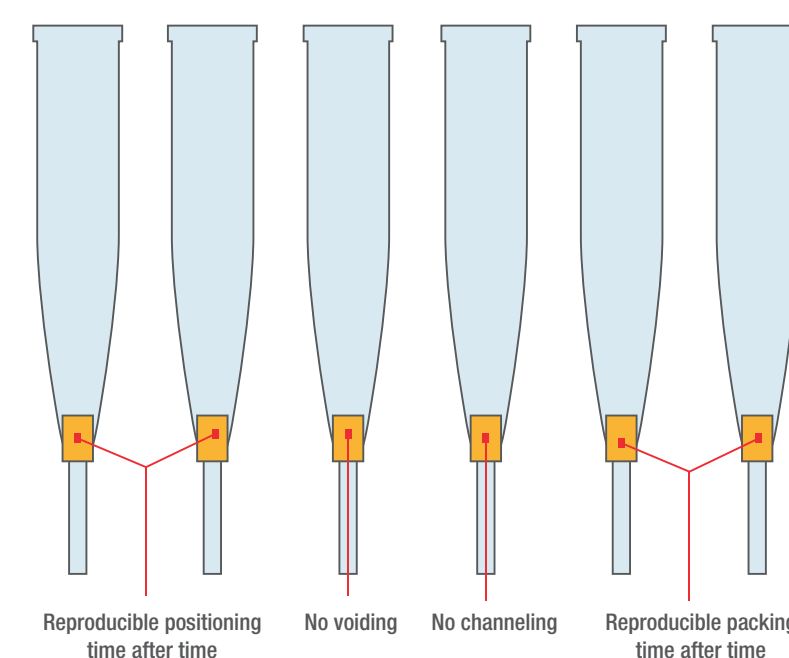
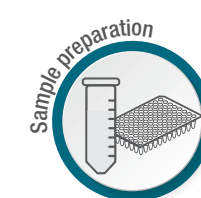
- Robustness – low analytical failure rates
- Ability to process low sample volumes
- High-sensitivity
- High-reproducibility
- Ease-of-use
- High-throughput processing
- Efficient and fast processes

Acidic compounds		Neutral compounds	Basic compounds	
Strong acids	Weak acids	Neutral compounds	Weak bases	Strong bases
(pK <sub>a</sub> < 2)	(pK <sub>a</sub> 2-4)		(pK <sub>a</sub> 8-10)	(pK <sub>a</sub> > 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™ SOLA $\mu$ ™ SPE range has the added benefit of being able to provide:

- Up to a 20 fold increase in sensitivity
- Ability to process samples restricted in volume
- Increased workflow efficiency and sample integrity



SOLA WAX	SOLA 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				
<b>Condition</b> 500 $\mu$ L methanol	<b>Condition</b> 500 $\mu$ L methanol	<b>Condition</b> 500 $\mu$ L methanol	<b>Condition</b> 500 $\mu$ L methanol	<b>Condition</b> 500 $\mu$ L methanol
<b>Equilibrate</b> 500 $\mu$ L water with 1% formic acid	<b>Equilibrate</b> 500 $\mu$ L water with 1% ammonium hydroxide	<b>Equilibrate</b> 500 $\mu$ L water	<b>Equilibrate</b> 500 $\mu$ L water with 1% formic acid	<b>Equilibrate</b> 500 $\mu$ L water with 1% ammonium hydroxide
<b>Load</b> 50-500 $\mu$ L of sample at 1 mL/min	<b>Load</b> 50-500 $\mu$ L of sample at 1 mL/min with 1% ammonium hydroxide	<b>Load</b> 50-500 $\mu$ L of sample at 1 mL/min	<b>Load</b> 50-500 $\mu$ L of sample at 1 mL/min with 1% formic acid	<b>Load</b> 50-500 $\mu$ L of sample at 1 mL/min with 1% ammonium hydroxide
<b>Wash 1</b> 500 $\mu$ L water	<b>Wash 1</b> 500 $\mu$ L water with 1% ammonium hydroxide	<b>Wash 1</b> 500 $\mu$ L 5% methanol in water	<b>Wash 1</b> 500 $\mu$ L water with 1% formic acid	<b>Wash 1</b> 500 $\mu$ L water with 1% ammonium hydroxide
<b>Wash 2</b> 500 $\mu$ L methanol	<b>Wash 2</b> 500 $\mu$ L methanol with 1% ammonium hydroxide		<b>Wash 2</b> 500 $\mu$ L methanol with 1% formic acid	<b>Wash 2</b> 500 $\mu$ L methanol with 1% ammonium hydroxide
<b>Elute</b> 200-500 $\mu$ L methanol with 1% ammonium hydroxide	<b>Elute</b> 200-500 $\mu$ L methanol with 1% formic acid	<b>Elute</b> 200-500 $\mu$ L methanol	<b>Elute</b> 200-500 $\mu$ L methanol with 1% ammonium hydroxide	<b>Elute</b> 200-500 $\mu$ L methanol with 1% formic acid

### SOLA $\mu$ SPE 96 Well Plates

Description	Bed weight	Well volume	Cat. no.	Quantity
SOLA $\mu$ HRP	2 mg	1 mL	60209-001	1 each
SOLA $\mu$ SCX	2 mg	1 mL	60209-002	1 each
SOLA $\mu$ SAX	2 mg	1 mL	60209-003	1 each
SOLA $\mu$ WCX	2 mg	1 mL	60209-004	1 each
SOLA $\mu$ 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

Description	Bed weight	Well volume	Cat. no.	Quantity
SOLA $\mu$ HRP	10 mg	1 mL	60109-001	100 pack
SOLA $\mu$ SCX	10 mg	1 mL	60109-002	100 pack
SOLA $\mu$ SAX	10 mg	1 mL	60109-003	100 pack
SOLA $\mu$ WCX	10 mg	1 mL	60109-004	100 pack
SOLA $\mu$ WAX	10 mg	1 mL	60109-005	100 pack

Find out more at [thermofisher.com/solaspe](https://thermofisher.com/solaspe)