

## Product Information

### Seppro® IgY 14 Tips

Catalog Number **S3824**

Storage Temperature 2–8 °C

## TECHNICAL BULLETIN

### Product Description

The Seppro® IgY 14 Tips are based on avian antibody (IgY)-antigen interactions and optimized buffers for sample loading, washing, eluting, and tip regeneration (available separately). The tips are specifically designed to remove fourteen highly abundant proteins from human biological fluids such as serum or plasma. The following proteins are depleted in a single step:

Albumin	IgG
$\alpha_1$ -Antitrypsin	IgA
IgM	Transferrin
Haptoglobin	$\alpha_2$ -Macroglobulin
Fibrinogen	Complement C3
$\alpha_1$ -Acid Glycoprotein (Orosomucoid)	
HDL (Apolipoproteins A-I and A-II)	
LDL (mainly Apolipoprotein B)	

The targeted highly abundant proteins are simultaneously removed by the immobilized specific IgYs when crude biological samples are passed through the tips. Selective immunodepletion of the highly abundant proteins provides enriched flow-through fractions of low abundance proteins for further study and downstream proteomics analysis. The removal of highly abundant proteins enables improved resolution and dynamic range for one-dimensional gel electrophoresis (1DGE), two-dimensional gel electrophoresis (2DGE), and liquid chromatography/mass spectrometry (LC/MS). The collected flow-through fractions may need to be concentrated dependent upon the downstream applications.

### Characteristics of IgY 14 Tips

The tips are designed for use with the automated Magtration® System SA-1 from Precision System Science USA. Twelve tips are simultaneously operated and twelve samples can be processed at once with no hand-on manipulations in ~65 minutes.

Tip Size: 0.5 ml bed volume

Tip capacity: 15  $\mu$ l of human serum/plasma  
(pooled normal human plasma or serum)

Operating temperature: 18–25 °C

Shipping Buffer: 1× Dilution Buffer with 0.02% sodium azide

Shipping temperature: 2–8 °C

Tip body materials: Polycarbonate tip and polyethylene frits

Usage: Tips may be used 30 times.

### Component

Seppro IgY 14 Tips 6 each  
(Catalog Number S3824)

### Reagents and Equipment Required for the Depletion Process, but Not Provided

10× Dilution Buffer Tris-Buffered Saline (TBS) - 100 mM Tris-HCl with 1.5 M NaCl, pH 7.4 (Catalog Number S4199)	1 × 200 ml
10× Neutralization Buffer 1 M Tris-HCl, pH 8.0 (Catalog Number S4449)	80 ml
Pre-filled Reagent Cartridges (Catalog Number S6574)	48 each
Corning® Spin-X® Centrifuge tube filters 0.45 $\mu$ m, pack of 100 (Catalog Number CLS8163)	1 × pack

### Precautions and Disclaimer

This product is for R&D use only, not for drug, household, or other uses. Please consult the Material Safety Data Sheet for information regarding hazards and safe handling practices.

### Storage/Stability

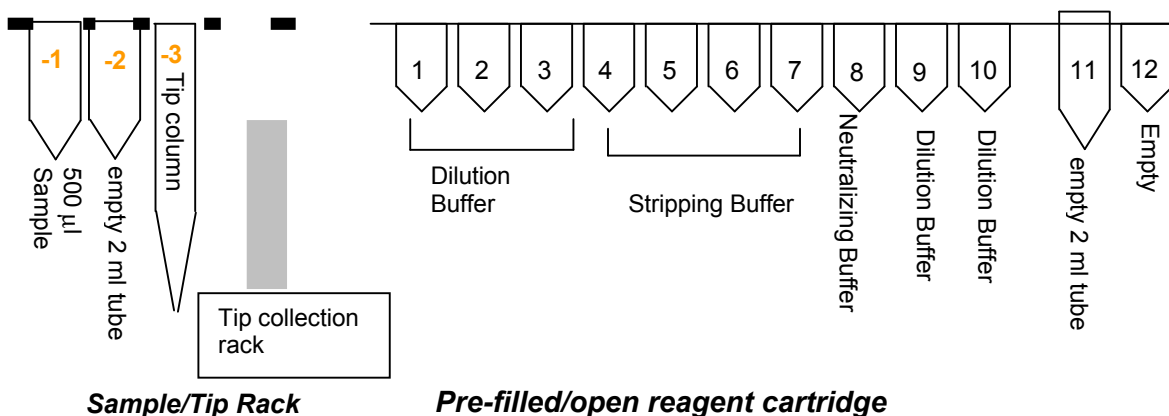
Store the tips at 2–8 °C. After use, equilibrate the tips with 1× Dilution Buffer containing 0.02% sodium azide and store them at 2–8 °C with the end-caps tightly sealed. **Do Not Freeze** the tips.

### Preparation Instructions

Preparation of 1× Dilution Buffer - Dilute 10× Dilution Buffer 10-fold with water. Use 1× Dilution Buffer for sample preparation.

Use 10× Neutralization Buffer to neutralize eluted bound proteins if analysis of bound proteins is desired.

Sample Preparation — It is not recommended to load an unfiltered serum or plasma directly onto the tip. Dilute the sample using 1× Dilution Buffer to a final volume of 500 µl. Samples may contain particulate materials, which can be removed by centrifugation at 10,000 rpm for 1 minute **or** with the use of a 0.45 µm spin filter with centrifugation for 1 minute at 9,000 × g. It is suggested to **avoid using reducing reagents, such as DTT, BME, or denaturing reagents, such as urea or guanidine-HCl** in the sample extracts.



The Buffer Arrangement in Pre-filled Cartridges and Tubes:

Position –1:	0.5 ml of unfractionated sample in 2 ml screw cap tube
Position –2:	Empty 2 ml screw cap tube
Position –3:	tip with tip holder
Wells 1–3:	1.0 ml of Dilution Buffer (Pre-filled Cartridge)
Wells 4–7:	1.0 ml of Stripping Buffer (Pre-filled Cartridge)
Well 8:	1.0 ml of Neutralizing Buffer (Pre-filled Cartridge)
Wells 9–10:	1.0 ml of Dilution Buffer (Pre-filled Cartridge)
Position 11:	Empty 2 ml screw cap tubes
Position 12:	Empty. Used for buffer retention.

**Note:** Buffers in wells are Pre-filled Cartridges.

### Procedure

**Notes:** Do not expose tips to organic solvents (like alcohols, acetonitrile, etc.), strong oxidizers, acids, or reducing agents and other protein denaturing agents.

Refer to Operation Manual for set-up of Magtration System SA-1 instrument, accessories, and disposables.

**Pretreatment of Tips** – “mock run” to remove any residual non-covalently bound IgY from the beads.

1. Turn on SA-1 and lift the door for reagent loading.

2. Load the Reagent Cartridges into the cartridge rack according to the number of the samples and tips to be used (SA-1 holds a maximum of 12 tips for 12 samples/per run).
3. Place the cartridge rack into the rack setting position on the stage.
4. Remove the top cap of the tip, leave the end cap with it, and put it at position –3 without tip sheath.
5. Set tip/tube rack into the instrument.
6. Insert the Tip Collection Rack at position –4

7. Close the door and follow the instructions on the display screen of the instrument to start the Mock Run (select 2) procedure for the process, see Table 1. It will take ~30 minutes to finish the process.

**Note:** After each run, remove both racks from the stage, replace all used cartridges and tubes. Take out tips with end caps on from tip collection rack and put them at position –3. Set racks back on the stage.

**Table 1.**

Mock Run

<Tip/tube rack>

Position –3	Seppro Tip in tip sheath
Position –2	Empty
Position –1	Empty

<Cartridge Rack>

Well 1	Equilibrium 1
Well 2	Equilibrium 2
Well 3	Equilibrium 3
Well 4	Elution 1
Well 5	Elution 2
Well 6	Elution 3
Well 7	Elution 4
Well 8	Neutralization
Well 9	Equilibrium 4
Well 10	Equilibrium 5
Position 11	Empty
Well 12	Retention liquids

Sample Depletion

- Put the tube containing the diluted human serum/plasma sample in a 2.0 screw cap tube at position –1. Put an empty 2 ml tube at position –2. Put a 2 ml tube at position 11.
- Start the depletion procedure (select 1) for sample processing (see Table 2).
- After completing the procedure, neutralize the eluted bound fractions by adding 100 µl of 10× Neutralization Buffer into the tube at position 11 (Elutions 1 & 2), well 4 (Elutions 3 & 4), and well 5 (Elutions 5 & 6); and 50 µl into well 6 (Elution 7) and well 7 (Elution 8).
- Transfer desired fractions (such as neutralized Elutions 3–8) to collection tubes. Store properly for downstream analysis.
- Store the tip with 1× Dilution Buffer at 2–8 °C. For long-term storage, use 1× Dilution Buffer with 0.02% sodium azide. **Do Not Freeze** the tips.

**Table 2.**

Depletion Run

<Tip/tube rack>

Position –3	Seppro Tip with end caps on
Position –2	Empty 2 ml tube. (After the process, this well contains ~0.5 ml of Wash 1).
Position –1	Sample 500 µl in 2 ml tube

<Cartridge Rack>

Well 1	Wash 2
Well 2	Wash 3
Well 3	Wash 4
Well 4	Elution 1 <b>(after the process, this well contains Elutions 3 and 4)</b>
Well 5	Elution 2 <b>(after the process, this well contains Elutions 5 and 6)</b>
Well 6	Elution 3 <b>(after the process, this well contains Elution 7)</b>
Well 7	Elution 4 <b>(after the process, this well contains Elution 8)</b>
Well 8	Neutralization
Well 9	Equilibrium 1, 2
Well 10	Equilibrium 3
Position 11	2 ml tube <b>(after the process, this well contains Elutions 1 and 2)</b>
Well 12	Retention liquids

Seppro is a registered trademark of Sigma-Aldrich® Biotechnology LP and Sigma-Aldrich Co.

Coming and Spin-X are registered trademarks of Corning, Inc.

Magtration is a registered trademark of Precision System Science Co., Ltd.

KR,DEC,MAM 12/08-1

Sigma brand products are sold through Sigma-Aldrich, Inc.

Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply. Please see reverse side of the invoice or packing slip.