

# Peptide/protein extraction protocol

Rev 1.0

When starting with serum or plasma samples, your antigen of interest may need to be extracted from contaminating peptides and proteins. To determine if extraction is required, spike the samples with a known amount of standard and check if they are accurately measured by the assay with and without extraction. An excess amount of antigen standard is provided to perform this procedure.

Extraction eliminates potentially interfering substances, such as bulk proteins and lipids. It can also be used to concentrate the sample so it falls within the measuring range. As with any purification technique, recovery of the desired substance is likely to be incomplete. Therefore, both optimization and quantification of the extraction procedure are recommended for accurate determinations.

## Required materials

- C18 Sep-Pak™ Column containing >50 mg of C18
- Binding Buffer ( 1% trifluoroacetic acid (TFA, HPLC grade))
- Elution Buffer (60% acetonitrile (HPLC grade), 1% TFA, and 39% distilled water)
- Protease inhibitor cocktail (e.g., Sigma P1860-1ML)

## Sample preparation guidelines

1. Collect plasma samples into EDTA collection tubes.
2. Centrifuge blood at  $1,600 \times g$  for 15 minutes at 4°C. Collect the top layer (plasma).
3. Add 0.5 µL of protease inhibitor cocktail for every 1 mL of plasma as soon as possible after collection.
4. Use the sample immediately or freeze in aliquots at -80°C or lower.

## Perform sample extraction (serum or plasma)

The following procedure is a generic protocol meant to help users with little experience in extracting samples. It is applicable to different biological fluids but cannot be thought of as an optimized protocol for any particular antigen.

1. Mix an equal amount of Binding Buffer to the plasma and vortex.
  - For **plasma** samples, mix an equal amount of Binding Buffer to the plasma and vortex.
  - For **tissue** samples, add 5mL of Binding Buffer for every gram of -80°C frozen tissue. Homogenize samples using a ultrasonic or Parr-type homogenizer.
  - For **CSF** samples, add add 1 mL of Binding Buffer for every 1 mL of CSF.
2. Centrifuge at  $6,000-17,000 \times g$  for 20 minutes at 4°C. Collect supernatant.
3. Equilibrate a C18 column by washing with 1 mL Elution Buffer.
4. Wash the C18 column with 3 mL of Binding Buffer three times.
5. Load the plasma/ Binding Buffer solution from step 2) onto the washed C18 column.
6. Apply light vacuum to the column (10 seconds/drop) and slowly wash the column with 3 mL of Binding Buffer twice. Discard the wash.
7. Apply light vacuum to the column (10 seconds/drop) and elute the peptide slowly with 3 mL of Elution Buffer. Collect eluate in a polypropylene tube.
8. Remove acetonitrile solvent in a centrifugal concentrator. Freeze-dry the resulting water/TFA solution to dryness.
9. Dissolve the residue in a suitable volume of 1X Assay Buffer provided with the kit.

## Important licensing information

This product may be covered by one or more Limited Use Label Licenses. By use of this product, you accept the terms and conditions of all applicable Limited Use Label Licenses.

## Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at [www.lifetechnologies.com/termsandconditions](http://www.lifetechnologies.com/termsandconditions). If you have any questions, please contact Life Technologies at [www.lifetechnologies.com/support](http://www.lifetechnologies.com/support)

## Product label explanation of symbols and warnings

	Catalog Number		Batch code		Temperature limitation		Use by		Manufacturer		Consult instructions for use		Caution, consult accompanying documents
--	----------------	---	------------	---	------------------------	---	--------	---	--------------	---	------------------------------	---	---

**DISCLAIMER:** LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

Corporate entity: Life Technologies | Carlsbad, CA 92008 USA | Toll Free in USA 1.800.955.6288

© 2017 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. Sep-Pak is a registered trademark of Waters Corporation.

For support visit [thermofisher.com/support](http://thermofisher.com/support) or contact [techsupport@lifetech.com](mailto:techsupport@lifetech.com)

[thermofisher.com](http://thermofisher.com)

2 January 2017