



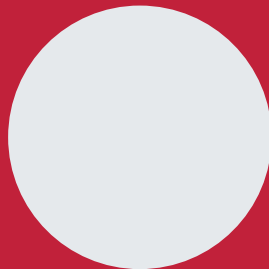
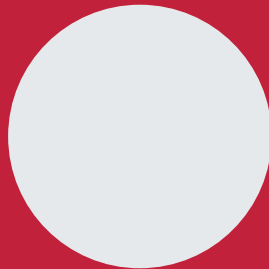
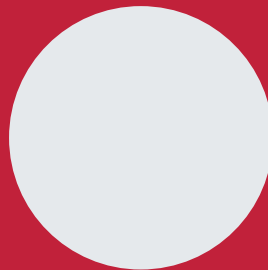
Innovating Epigenetic Solutions

# CaptEV reagent

**CaptEV: EV pre-enrichment reagent**

Cat. No. **C28030001** (cell culture)

Cat. No. **C28030002** (serum/plasma)



the 1990s, the number of people in the UK who are employed in the public sector has increased from 10.5 million to 13.5 million, and the number of people in the public sector who are employed in health care has increased from 2.5 million to 3.5 million (Department of Health 2000).

There are a number of reasons for the increase in the number of people employed in the public sector. One reason is that the public sector has become a major employer in the UK. Another reason is that the public sector has become a major employer in the health care sector. A third reason is that the public sector has become a major employer in the social care sector. A fourth reason is that the public sector has become a major employer in the education sector.

The increase in the number of people employed in the public sector has led to a number of changes in the way that the public sector is organized. One change is that the public sector has become more decentralized. Another change is that the public sector has become more market-oriented. A third change is that the public sector has become more customer-oriented. A fourth change is that the public sector has become more performance-oriented.

The changes in the way that the public sector is organized have led to a number of challenges for the public sector. One challenge is that the public sector has become more complex. Another challenge is that the public sector has become more competitive. A third challenge is that the public sector has become more demanding. A fourth challenge is that the public sector has become more demanding.

The challenges that the public sector faces are a result of the changes in the way that the public sector is organized. The challenges that the public sector faces are a result of the changes in the way that the public sector is organized. The challenges that the public sector faces are a result of the changes in the way that the public sector is organized. The challenges that the public sector faces are a result of the changes in the way that the public sector is organized.

The challenges that the public sector faces are a result of the changes in the way that the public sector is organized. The challenges that the public sector faces are a result of the changes in the way that the public sector is organized. The challenges that the public sector faces are a result of the changes in the way that the public sector is organized. The challenges that the public sector faces are a result of the changes in the way that the public sector is organized.

The challenges that the public sector faces are a result of the changes in the way that the public sector is organized. The challenges that the public sector faces are a result of the changes in the way that the public sector is organized. The challenges that the public sector faces are a result of the changes in the way that the public sector is organized. The challenges that the public sector faces are a result of the changes in the way that the public sector is organized.

The challenges that the public sector faces are a result of the changes in the way that the public sector is organized. The challenges that the public sector faces are a result of the changes in the way that the public sector is organized. The challenges that the public sector faces are a result of the changes in the way that the public sector is organized. The challenges that the public sector faces are a result of the changes in the way that the public sector is organized.

# Content

Introduction

Required materials not provided

Remarks before starting

Protocol

Related products

## Introduction

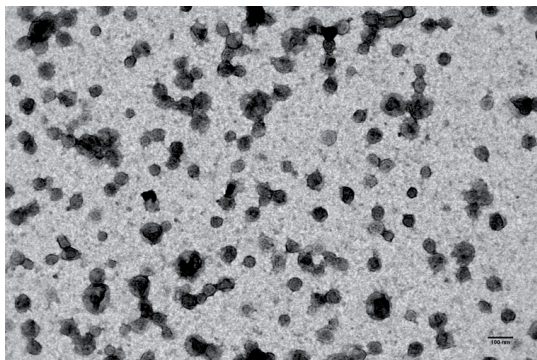
CaptEV precipitation reagent has been specifically designed as a pre-enrichment tool prior to extracellular vesicle purification. This reagent is ideally suited for total extracellular vesicle (EV) recovery for one biofluid that can later be processed by immunocapture using Diagenode's ExoIP™ kit.

The CaptEV reagent has been designed to precipitate all vesicles within a single sample to avoid bias. After precipitation, the exosome-rich suspension can be further purified by several means. The reagent also contains chemicals that support downstream immunocapture by facilitating exosome-bead complex formation. These chemicals are simple to wash away after the exosome isolation providing a highly enriched sample. As shown in Figure 1, exosomes are not trapped in sticky polymer and are efficiently isolated due to the CaptEV.

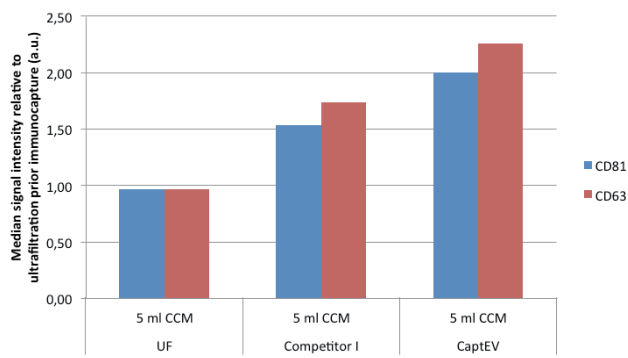
The CaptEV reagent is straightforward to use and allows for reproducibility between experiments.

### Advantages of CaptEV:

1. Full recovery of EVs
2. Enhanced immunocapture
3. Better signal coverage in FACS than with other reagents
4. Ease of use
5. Reproducibility
6. Intact and preserved exosomes (not sticky)



**Figure 1:** TEM picture of an exosome preparation obtained with Diagenode ExoIP anti-CD81 kit with elution of beads and Diagenode CaptEV reagent as a pre-enrichment tool



**Graph 1:** Median signal intensity from the flow cytometry analysis of the exosome capture efficiency corresponding to different pre-enrichment ways prior immunocapture (ExoIP™)

## Required materials not provided

- Phosphate Buffered Saline (PBS)
- Pipetous
- Serological pipettes
- Micropipettes and corresponding tips: 20µl – 200µl – 1000µl
- 15ml and/or 50ml tubes for cell culture supernatant collection
- Centrifuge
- Gloves
- Crushed ice

## Remarks before starting

CaptEV reagent for cell culture supernatant or for serum and plasma has been designed to be used as a pre-enrichment tool prior to EV purification.

It has been specially designed to work with Diagenode's ExoIP™ immunocapture kits. However, other purification may be used if desired.

Moreover, concerning the cell culture supernatant; this reagent works well with conditioned cell culture supernatant. In normal conditions of use, exosome-depleted FBS complemented cell culture media can be used as starting material.

The use of CaptEV reagent for cell culture supernatant has only been validated for use with samples sourced from cell culture supernatant. Other sample types have not been tested.

## Protocol

The reagent is supplied in a 2X concentrated form. One (1) volume of reagent has to be mixed with one (1) volume of biofluid.

1. Harvest biofluid containing the extracellular vesicles of interest.
2. For cell culture supernatant, spin it at 3000g for 30 minutes at +4°C in order to eliminate floating cells and cells debris.
3. Filter biofluid gently through a 0.2 µm cut-off to eliminate large vesicles such as apoptotic bodies and most of the microvesicles.
4. Mix one volume of cleared, conditioned cell culture supernatant or serum and plasma with one volume of precipitation reagent.

The precipitation reaction can be easily scaled up or down depending on the needs of the final user. The precipitation reagent is more efficient than an ultrafiltration column when used in combination with immunocapture (Diagenode ExoIP™ kit). Thus, a protocol set up previously for concentrating cell culture supernatant can be used as is for precipitation before immunocapture.

5. Mix the precipitation mixture by gently inverting the reaction tube several times.
6. Incubate 1 hour at room temperature (RT) or at +4°C overnight (16 hours) under mild agitation.
7. Spin the precipitation mixture at 1500g for 30 minutes at +4°C.
8. Discard the supernatant.
9. Resuspend the pellet with the desired volume of PBS before downstream processing.

Typically, the pellet should be very easily resuspendable except if the cell culture supernatant contains poorly depleted serum.

**Notice:** If an exosome-rich pellet is processed by immunocapture (ExoIP™) afterwards, it is advised to keep the resuspension volume under 500 µl of PBS. Too large of a volume (→ 1 ml) is not beneficial, since bead-binding kinetics is dependent on the reaction volume.

10. Process the exosome-enriched suspension into immunocapture with the ExoIP™ kit or store it at -80°C if not used immediately.

At this stage, the exosome preparation is far from being pure. Further purification such as immunocapture (ExoIP™) of the target or gel filtration (e.g. Diagenode EVCleaner) need to be applied. Moreover, the exosome-rich preparation obtained by precipitation contains chemicals that enhance downstream immunocapture but may not be suitable for

direct processing by other techniques that require a very pure sample.

## Related products

| Product name | Cat. No.            |
|--------------|---------------------|
| C28030001    | CaptEV cell culture |
| C28030002    | CaptEV serum/plasma |
| C28020001    | EVCleaner           |

## **FOR RESEARCH USE ONLY.**

**Not intended for any animal or human therapeutic or diagnostic use.**

© 2016 Diagenode SA. All rights reserved. No part of this publication may be reproduced, transmitted, transcribed, stored in retrieval systems, or translated into any language or computer language, in any form or by any means: electronic, mechanical, magnetic, optical, chemical, manual, or otherwise, without prior written permission from Diagenode SA (hereinafter, "Diagenode"). The information in this guide is subject to change without notice. Diagenode and/or its affiliates reserve the right to change products and services at any time to incorporate the latest technological developments. Although this guide has been prepared with every precaution to ensure accuracy, Diagenode and/or its affiliates assume no liability for any errors or omissions, nor for any damages resulting from the application or use of this information. Diagenode welcomes customer input on corrections and suggestions for improvement.

# **NOTICE TO PURCHASER**

## **LIMITED LICENSE**

The information provided herein is owned by Diagenode and/or its affiliates. Subject to the terms and conditions that govern your use of such products and information, Diagenode and/or its affiliates grant you a nonexclusive, nontransferable, non-sublicensable license to use such products and information only in accordance with the manuals and written instructions provided by Diagenode and/or its affiliates. You understand and agree that except as expressly set forth in the terms and conditions governing your use of such products, that no right or license to any patent or other intellectual property owned or licensable by Diagenode and/or its affiliates is conveyed or implied by providing these products. In particular, no right or license is conveyed or implied to use these products in combination with any product not provided or licensed to you by Diagenode and/or its affiliates for such use.

Limited Use Label License: Research Use Only The purchase of this product conveys to the purchaser the limited, non-transferable right to use the product only to perform internal research for the sole benefit of the purchaser. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. This product is for internal research purposes only and is not for use in commercial applications of any kind, including, without limitation, quality control and commercial services such as reporting the results of purchaser's activities for a fee or other form of consideration. For information on obtaining additional rights, please contact [info@diagenode.com](mailto:info@diagenode.com).

## **TRADEMARKS**

The trademarks mentioned herein are the property of Diagenode or their respective owners. Bioanalyzer is a trademark of Agilent Technologies, Inc. Agencourt and AMPure are registered trademarks of Beckman Coulter, Inc. Illumina is a registered trademark of Illumina Inc; Qubit is a registered trademark of Life Technologies Corporation.



Innovating Epigenetic Solutions

**DIAGENODE  
HEADQUARTERS**

**DIAGENODE S.A.  
BELGIUM | EUROPE**

LIEGE SCIENCE PARK  
Rue Bois Saint-Jean, 3  
4102 Seraing - Belgium  
Tel: +32 4 364 20 50  
Fax: +32 4 364 20 51  
orders@diagenode.com  
info@diagenode.com

**DIAGENODE INC.  
USA | NORTH AMERICA**

400 Morris Avenue, Suite #101  
Denville, NJ 07834  
Tel: +1 862 209-4680  
Fax: +1 862 209-4681  
orders.na@diagenode.com  
info.na@diagenode.com

**FOR A COMPLETE LISTING OF  
DIAGENODE'S INTERNATIONAL  
DISTRIBUTORS VISIT:**

[http://www.diagenode.com/company/  
distributors.php](http://www.diagenode.com/company/distributors.php)  
For rest of the world, please contact  
Diagenode sa.  
[www.diagenode.com](http://www.diagenode.com)