

# EXO GAG

## CEREBRO SPINAL FLUID

Next generation of Exosome isolation  
for biomarker discovery in liquid biopsy

### EXOSOME PURIFICATION KIT FOR CEREBRO SPINAL FLUID

#### USER GUIDE

#### STORAGE

---

All components can be stored at room temperature.

#### PRODUCT COMPONENT

---

EXO GAG - CSF™ isolation kit 20ml (40 samples). 1 x User guide.

Not supplied: 1.5ml microcentrifuge collection tubes.

**EXO GAG - CSF is a specific, quick and inexpensive method to optimize the process of exosome isolation.**

#### PRODUCT INFORMATION

---

The EXO GAG - CSF precipitation reaction is based on the interaction between the precipitation solution and glycosaminoglycans located on exosomes surface. After a simple incubation, the exosomes can be isolated by a short centrifugation.

#### WHAT IS EXO GAG - CSF

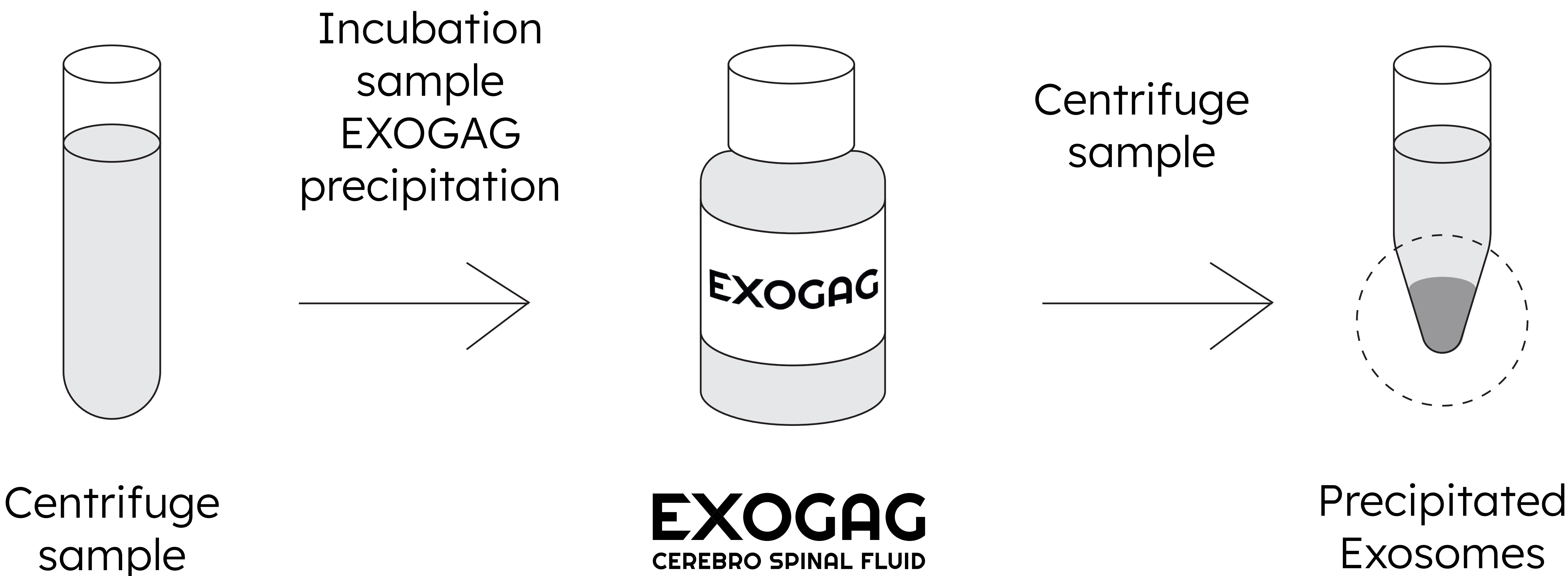
---

EXO GAG - CSF technology allows the isolation of exosomes from cerebro spinal fluid based on the affinity of exosomes to the precipitation reagent, which allows their isolation from a complex sample. EXO GAG - CSF is a patented exosome purification method that allows the isolation of exosomes from a cerebro spinal fluid sample with a minimal amount of co-precipitated material, such as protein or genetic material (DNA, RNA or microRNA) which makes it an ideal product for biomarkers research and their transfer to the clinic.



|                                  |      |             |                     |                              |
|----------------------------------|------|-------------|---------------------|------------------------------|
| High sensibility and specificity | Fast | Inexpensive | Small sample needed | No specific equipment needed |
|----------------------------------|------|-------------|---------------------|------------------------------|

|                      |              |                   |
|----------------------|--------------|-------------------|
| SAMPLE               | SAMPLE, vol. | EXOAG - CSF, vol. |
| Cerebro Spinal Fluid | 250µl        | 500µl             |



## PROTOCOL

### EXOAG - CSF Exosomes Isolation Protocol.

- 1. Collect CSF sample.** Samples can be frozen until the moment they are used; if the samples have been frozen, thaw and temper them before processing.
- 2. Centrifuge the sample** at 2000 x g for 5' to remove cells and cell debris.
- 3. Transfer the supernatant** to a new tube and discard the pellet of possible cell debris.
- 4. Add the volume of sample** to isolate exosomes to a new tube and add twice the volume of EXOGAG - CSF precipitation reagent, as shown in the table.
- 5. Mix the sample** and EXOGAG - CSF precipitation reagent by inverting the tube or vortexing to homogenize the final solution (the solution will have a characteristic blue colour).
- 6. Incubate the sample** for 5' at 4°C.
- 7. Centrifuge the sample** at 3000 x g; 15' at 4°C.
- 8. Remove the supernatant** being careful not to remove the pellet containing the exosomes (this pellet will be dark blue).
- 9. Resuspend the exosomes** in the appropriate buffer (pipetting repeatedly up and down), depending on the technique.