

## ab286884 – Lyophilized Exosome Standard (100 µg, Human Urine)

For use as a standard positive control for immunocapture performance evaluation.

For research use only - not intended for diagnostic use.

For overview, typical data and additional information please visit:

<http://www.abcam.com/ab286884>

### Storage and Stability

On receipt product should be stored at +4°C. Upon opening, use kit within 36 months, provided lyophilized exosome standard has not been reconstituted.

### Materials Supplied

Item	Quantity 2 vials	Quantity 4 vials	Quantity 6 vials	Storage Condition
Lyophilized Exosome Standard	2 x 100 µg	4 x 100 µg	6 x 100 µg	+4 °C

### Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully utilize this assay:

- Deionized water

### Reagent Preparation

Lyophilized Exosome Standard: Reconstitute lyophilized exosome standard by adding deionized water. Reconstituted exosome standards are not suitable for long term conservation at room temperature, use them within 2 hours after reconstitution. The remaining reconstituted standard stock solution should be aliquoted into polypropylene vials (preferably low binding) and stored at -20°C for up to one month or at -80°C for up to six months. Strictly avoid repeated freeze-and-thaw cycles.

### Assay Protocol

**Δ Note:** Different volumes of deionized water for exosomes reconstitution can be chosen by the users in according with the desired final concentration.

1. Reconstitute lyophilized exosome standard by adding deionized water. 100 µl for lyophilized Standard 100 µg, to get a final concentration of 1 µg/µl.
2. Suspend exosomes pipetting the solution up and down 10-15 times, avoiding bubbles.
3. Vortex the reconstituted standard for 60 secs.
4. Briefly centrifuge the tubes containing the standard to ensure that the solution is collected at the bottom of the tube.
5. Pipette the solution up and down 10 times, avoiding the introduction of bubbles. After this step, the standard is ready to use.

### **Technical Support**

Copyright © 2021 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

For all technical or commercial enquiries please go to:

[www.abcam.com/contactus](http://www.abcam.com/contactus)

[www.abcam.cn/contactus](http://www.abcam.cn/contactus) (China)

[www.abcam.co.jp/contactus](http://www.abcam.co.jp/contactus) (Japan)