

# Examining Exosomes on Simple Western

## Introduction

Exosomes are small endosome-derived lipid nanoparticles approximately 30-200 nm in diameter that are actively secreted by exocytosis by most living cells<sup>1</sup>. Exosome release can occur constitutively or upon induction, under pathological or non-pathological conditions, and in a dynamic, regulated and functionally-relevant manner. The molecular composition and quantity of released exosomes depend on the state of the parent cell. Exosomes are isolated from diverse cell lines and primary cultures, as well as from biological fluids like serum, plasma and saliva. Given this diversity, exosomes have pleiotropic physiological functions, and it is becoming increasingly clear that they have an important role in many pathological conditions such as cancer, infectious diseases, and neurodegenerative diseases. Hence, exosomes are a prime target in liquid biopsy methods<sup>2</sup>. Exosomes may also be engineered as promising vesicles for drug delivery<sup>3</sup>.

Due to their tiny size, isolating exosomes from complex biological matrices can be a challenge, and yields are often low. Here, Simple Western™ is an ideal solution for studying exosomes because it requires only 3 µL of sample and offers picogram-level sensitivity, reducing the number of exosomes typically required for traditional Western blot. Characterization of exosomes can subsequently be performed using appropriate detection antibodies against exosome-associated antigens. These target antigens may be generic exosome markers or specific to the cell or fluid sample from which they originate. In this protocol, we use Simple Western to examine both generic and sample-dependent targets using commercially-available exosome standards and antibodies.

## Materials

### ProteinSimple

- Instrument: Peggy Sue, Sally Sue, Jess or Wes
- 12–230 kDa Separation Module for Peggy Sue or Sally Sue (PN [SM-S001](#))
- 12–230 kDa Separation Module for Jess or Wes (PN [SM-W004](#))
- Anti-Rabbit Detection Module for Jess, Wes, Peggy Sue or Sally Sue (PN [DM-001](#))
- RIPA Lysis Buffer (PN [040-483](#))

### Exosome Standards

SOURCE	PART NUMBER	SOURCE	PART NUMBER
Plasma (Human)	NBP2-49838	DAUDI Cells (Human)	NBP2-49860
Serum (Human)	NBP2-49827	HCT116 Cells (Human)	NBP2-49854
<b>A549 Cells (Human)</b>	<b>NBP2-49862</b>	K562 Cells (Human)	NBP2-49864
Saliva (Human)	NBP2-49842	MM1 Cells (Human)	NBP2-49847
Urine (Human)	NBP2-49840	PC3 Cells (Human)	NBP2-49856
BLCL21 Cells (Human)	NBP2-49849	SK-N-SH Cells (Human)	NBP2-49852
BPH-1 Cells (Human)	NBP2-49858	U87-MG Cells (Human)	NBP2-49844
COLO1 Cells (Human)	NBP2-49845	B16F10 Cells (Mouse)	NBP2-49866

**TABLE 1.** Exosome standards available from [Novus Biologicals](#). Highlighted in bold are the standards used in this study.

### Antibodies

TARGET	VENDOR	PART NUMBER	DILUTION USED IN THIS STUDY
Alix	Novus Biologicals	NBP1-49701	1:10
Annexin A2	Cell Signaling Technologies	8235	1:250
Annexin V	Novus Biologicals	NB100-1930	1:250
CD63	Abcam	ab68418	1:10
CD9	Cell Signaling Technologies	D801A	1:10
Flotillin-1	Novus Biologicals	NBP1-79022	1:250
HER1/EGFR	Cell Signaling Technologies	2646	1:50
HSPA8/HSC70	Cell Signaling Technologies	8444	1:50
TSG101	Sigma	T5826	1:10

**TABLE 2.** Antibodies against exosome targets validated on Simple Western. All antibodies are sourced from rabbit. These targets were chosen from the list of top exosome markers from [ExoCarta](#).

## Methods

1. For analysis on Simple Western, lyse the exosome standards directly in 1X RIPA on ice for 15 minutes and then proceed with assay setup. Typically, a final exosome-derived lysate sample of 0.2 mg/mL was prepared per the product insert of the Separation Module.

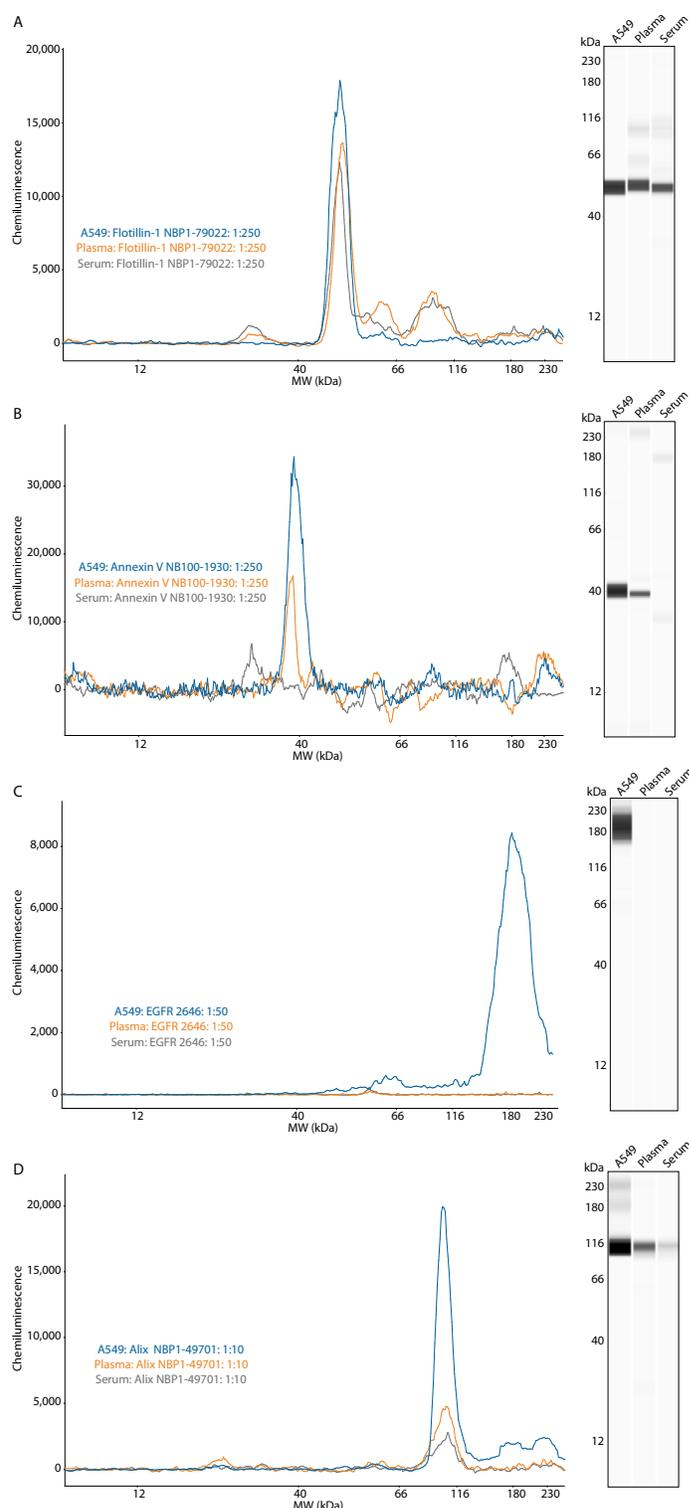
If you need exosomes for additional applications, first resuspend the exosomes in deionized water per the instructions provided in the product insert of the exosome standards prior to lysis in RIPA buffer.

Additional denaturation conditions should be considered when working with specific classes of proteins. For example, when working with membrane proteins, follow the guidelines in [Analyzing Integral Membrane Proteins with Simple Western](#). This protocol also describes treatment of glycosylated proteins with PNGase F.

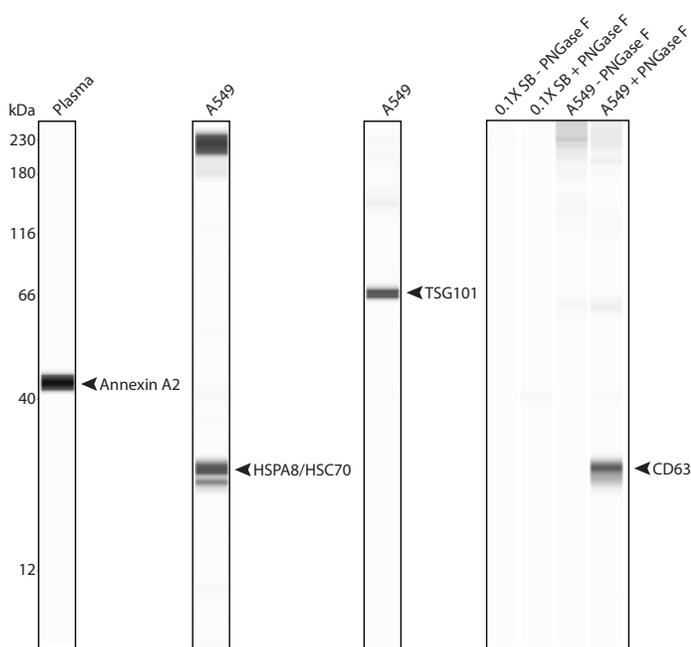
2. Dilute the antibodies in Antibody Diluent 2.
3. Follow the default sample preparation and assay conditions for Peggy Sue, Sally Sue, Jess or Wes.

## Results

We examined three exosome standards from Novus Biologicals (**Table 1**) derived from human plasma, human serum and A549 cells, which are adenocarcinomic human alveolar basal epithelial cells. Using these standards, we validated antibodies to common exosome targets listed in **Table 2**. Certain targets, such as the integral membrane protein Flotillin-1, are canonical exosome markers and ubiquitously expressed in most exosomes. As expected, Flotillin-1 was found in all exosome standards we tested (**Figure 1A**). Conversely, some proteins are specific to the origin of the exosome. For example, Annexin V, found in the plasma membrane, is a marker of apoptosis and anticoagulant protein in the blood. As such, it was found in A549 and plasma exosomes, but not found in serum-derived exosomes (**Figure 1B**). Likewise, EGFR is specific for A549-derived exosomes and it was not detected in exosomes derived from serum or plasma (**Figure 1C**). Finally, CD63 (Alix) was detected in all three exosome standard preparations, but had higher overall expression in A549-derived exosomes (**Figure 1D**).



**FIGURE 1.** Detection of exosome targets by Simple Western. (A) Flotillin-1 (B) Annexin V (C) EGFR and (D) Alix. Electropherograms are shown (left), as well as lane view (right).



**FIGURE 2.** Detection of additional exosome targets by Simple Western. Shown here is the lane view of each target detected on exosome standards from plasma and A549 cells. CD63 gave a signal only upon treatment with PNGase F. The lanes labeled with 0.1X SB contain sample buffer but lack exosome standard. For more information on PNGase F treatment, refer to [Analyzing Integral Membrane Proteins with Simple Western](#).

The other antibodies to common exosome targets listed in **Table 2** were validated on Simple Western (**Figure 2**). Interestingly, a signal for CD63 appeared only upon treatment of the exosome standard with the deglycosylation enzyme PNGase F (**Figure 2**), which highlights the importance of taking into consideration the specific class of a target protein, for example whether it is glycosylated or membrane-associated.

## Conclusion

When reconstituted to a concentration of 1 mg/mL, the exosome standards in **Table 1** have at least  $1 \times 10^{10}$  particles/mL. Therefore, at a final sample concentration of 0.2 mg/mL, the theoretical count of exosomes analyzed per well with a 3  $\mu$ L loading volume is approximately 6 million. However, this is certainly not the lower limit of detection; we were able to measure a robust signal of Flotillin-1 with as little as 0.03 mg/mL (900,000 particles) of pure exosome sample. Using these standards, we validated antibodies to common exosome targets on Simple Western (**Table 2**). These results show that Simple Western is an ideal method for analysis of exosomes where sample volume may be of critical importance.

## References

1. Exosomes, D M Pegtel and S J Gould, *Annual Review of Biochemistry*, 2019, 88:487–514.
2. Extracellular vesicles – biomarkers and effectors of the cellular interactome in cancer, J Rak, *Frontiers in Pharmacology*, 2013, 4:21.
3. Exosomes as reconfigurable therapeutic systems, R Conlan, S Pisano, M Oliveira, M Ferrari and I Mendes Pinto, *Trends in Molecular Medicine*, 2017, 23(7):636–650.