

KEY FINDING Preformed loadable SM-102 lipid nanoparticles enabled effective delivery of nucleic acid cargo to ovarian cancer and macrophage cell lines to modulate ENPP1 activity and STING activation.

INTRODUCTION

Ovarian cancer is a complex and heterogeneous disease with various subtypes and diverse molecular mechanisms of pathogenesis. Among different forms, epithelial ovarian cancer is the most common, representing up to 90% of total reported cases. Recent studies have shown that ENPP1 may play a role in cancer cell proliferation, migration, and invasion in ovarian and other types of cancer. ENPP1 is a glycosylated, type II transmembrane protein acting as a pyrophosphatase and phosphodiesterase with broad specificity. ENPP1 substrates include cyclic dinucleotides such as 2'3'-cGAMP, which binds to and activates STING, leading to the activation of the STING-dependent type 1 interferon (IFN) pathway and stimulation of antitumor activity.²⁻⁴ ENPP1 is frequently upregulated during tumorigenesis where its cGAMP hydrolytic activity contributes to an immunosuppressive tumor microenvironment.⁵ In this study we used preformed lipid nanoparticles (LNPs) to modulate ENPP1 activity and STING activation in ovarian cancer cell lines and THP-1 macrophages, respectively.

Objectives

- Knockdown ENPP1 in PA-1, Caov-3, and SKOV3 ovarian cancer cell lines via loadable SM-102 LNPs.
- Assess whether reduced ENPP1 activity leads to an increase in 2'3'-cGAMP production.
- Test whether ENPP1 is present on extracellular vesicles (EVs) in ovarian cancer cell lines.
- Demonstrate utility of loadable, lyophilized SM-102 LNPs as a research tool in delivering nucleic acid cargo to different cell types.

MATERIALS & METHODS

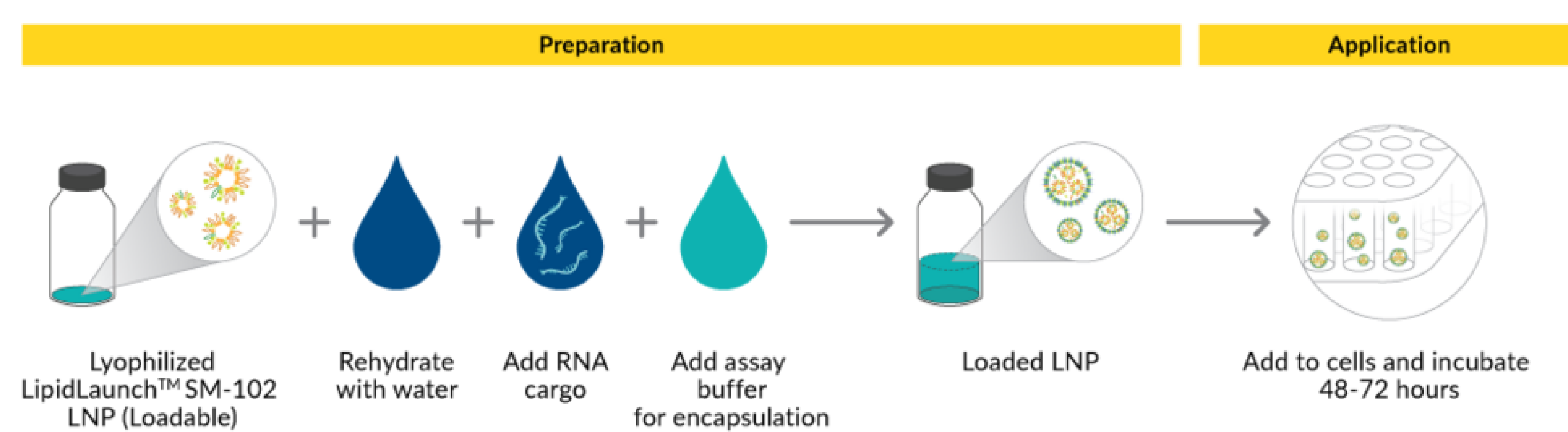


FIGURE 1 – Overview of cargo loading and transfection via loadable LNPs.

PA-1, Caov-3, and SKOV3 cells were seeded into 96- or 24-well plates at densities of 4,000 or 20,000 cells per well, respectively, 24 hrs prior to treatments. ENPP1 siRNA (50 to 250 ng per well) was loaded into preformed SM-102 LNPs, and cells were transfected for up to 72 hrs prior to assay.

ENPP1 activity was determined using Cayman's ENPP1/ENPP3 Cell-Based Activity Assay Kit (Item No. 702080). For 2'3'-cGAMP quantitation, PA-1 cells were refreshed with serum-free media 72 hrs after siRNA treatment and incubated for 24 hrs prior to detection of 2'3'-cGAMP in the conditioned media using Cayman's 2'3'-cGAMP ELISA Kit (Item No. 501700). EV fractions were prepared from conditioned media using Cayman's Extracellular Vesicle Isolation Kit for tissue cultures (Item No. 702630).

MATERIALS & METHODS CONTINUED

2'3'-cGAMP (Item No. 19887) or Cy5-labeled 2'3'-cGAMP (Biolog) was delivered to THP-1 macrophages via preformed loadable LNPs. Activation of STING was assessed after 24 hrs via immunoblotting (STING mAb; Item No. 17856). Uptake of labeled 2'3'-cGAMP was determined via fluorescence imaging (Cytation™ 5, Agilent). IL-1β and IL-6 were quantified in THP-1 conditioned media using Cayman's Interleukin-1β (human) ELISA Kit (Item No. 583311) and Interleukin-6 (human) ELISA Kit (Item No. 501030).

RESULTS

ENPP1 mRNA levels, protein levels, and enzymatic activity are suppressed in PA-1, SKOV3, and Caov-3 cells following loadable LNP-mediated siRNA knockdown

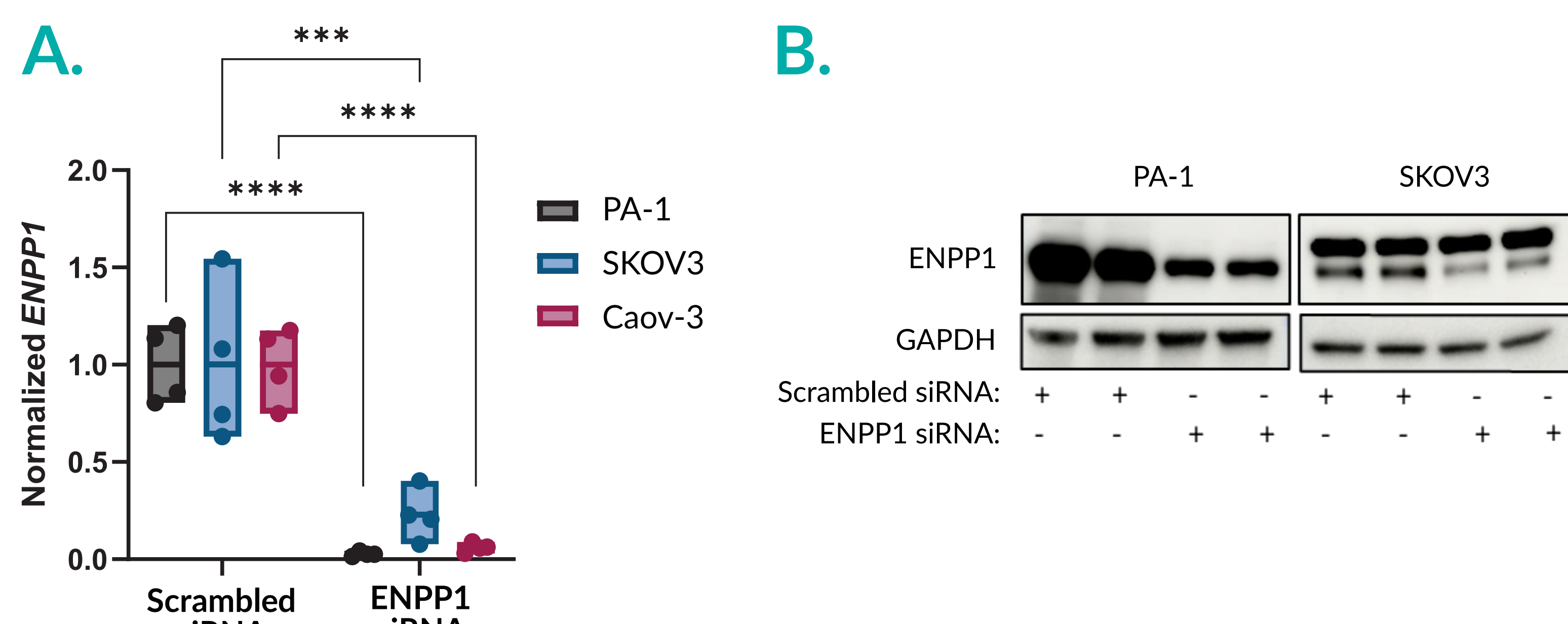


FIGURE 2 – Expression levels of ENPP1 in various ovarian cancer cell lines following siRNA-mediated knockdown.

The mRNA levels of ENPP1 in PA-1, SKOV3, and Caov-3 cell lines is significantly reduced following transfection with siRNA-loaded vs. scrambled siRNA-loaded SM-102 LNPs (A). Protein expression was determined via Western blotting (B). N=4, ****<0.0001.

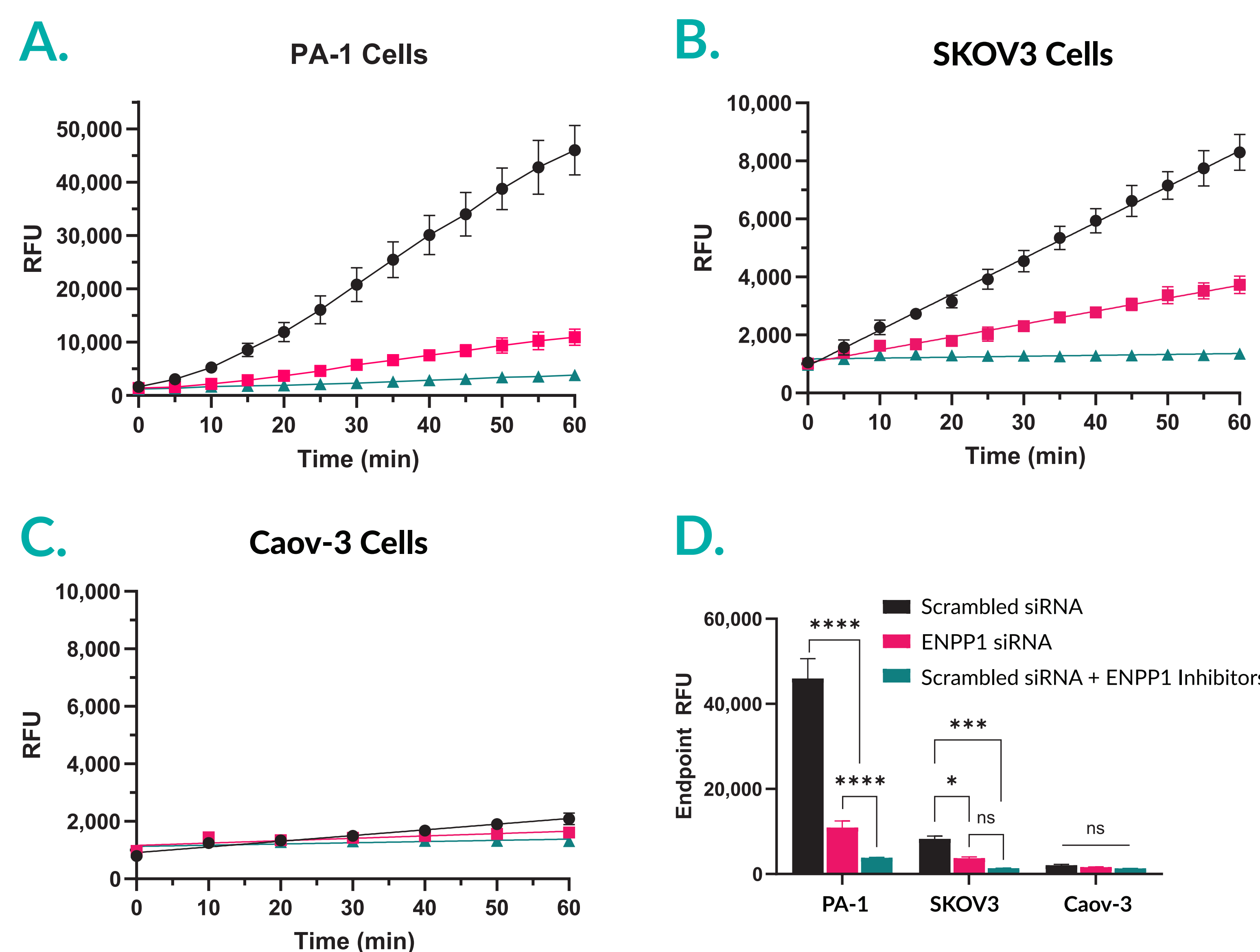


FIGURE 3 – ENPP1 activity in various ovarian cancer cell lines following siRNA-mediated knockdown.

ENPP1 activities were observed over a period of 60 minutes for PA-1 (A), SKOV3 (B), and Caov-3 (C). The endpoints after the 1-hr incubation are shown in (D). N=5, ****p<0.0001, ***p<0.001, *p<0.05.

RESULTS CONTINUED

Extracellular 2'3'-cGAMP levels are increased in ENPP1-depleted PA-1 cells

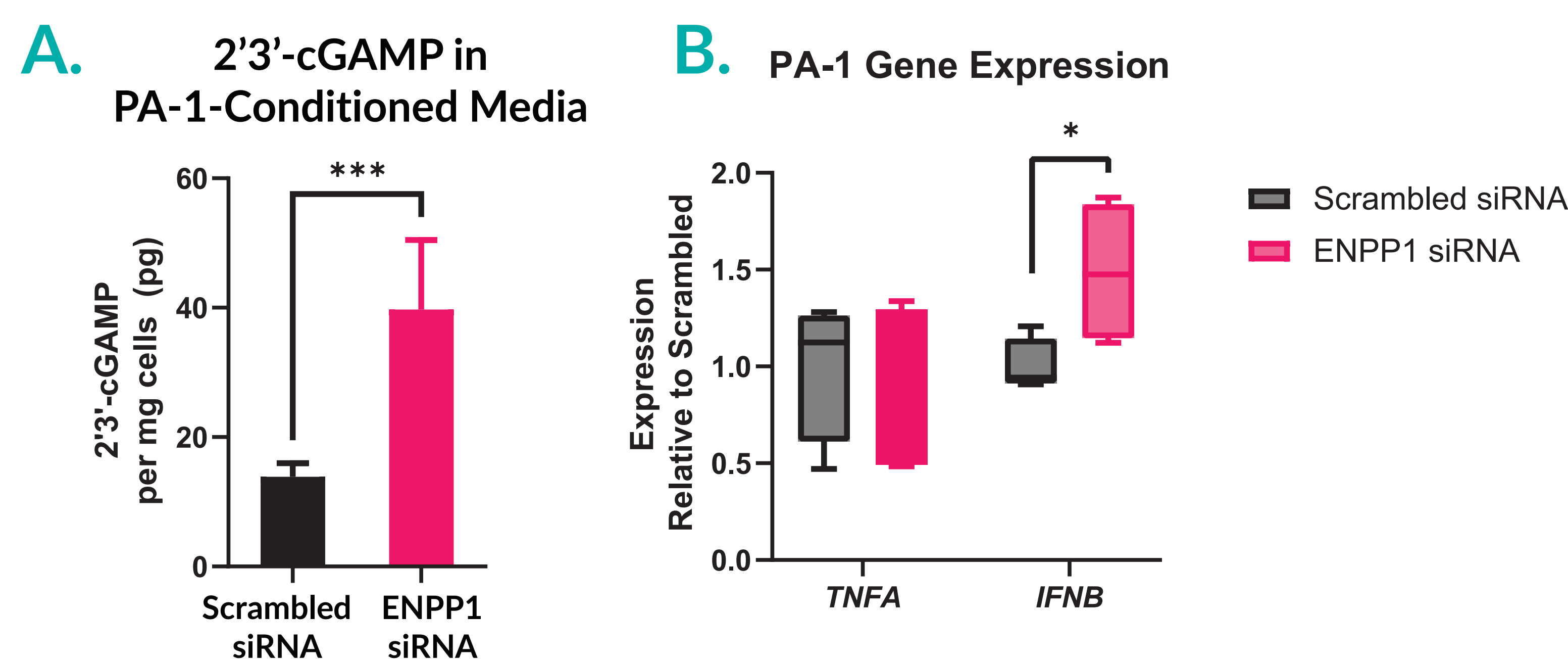
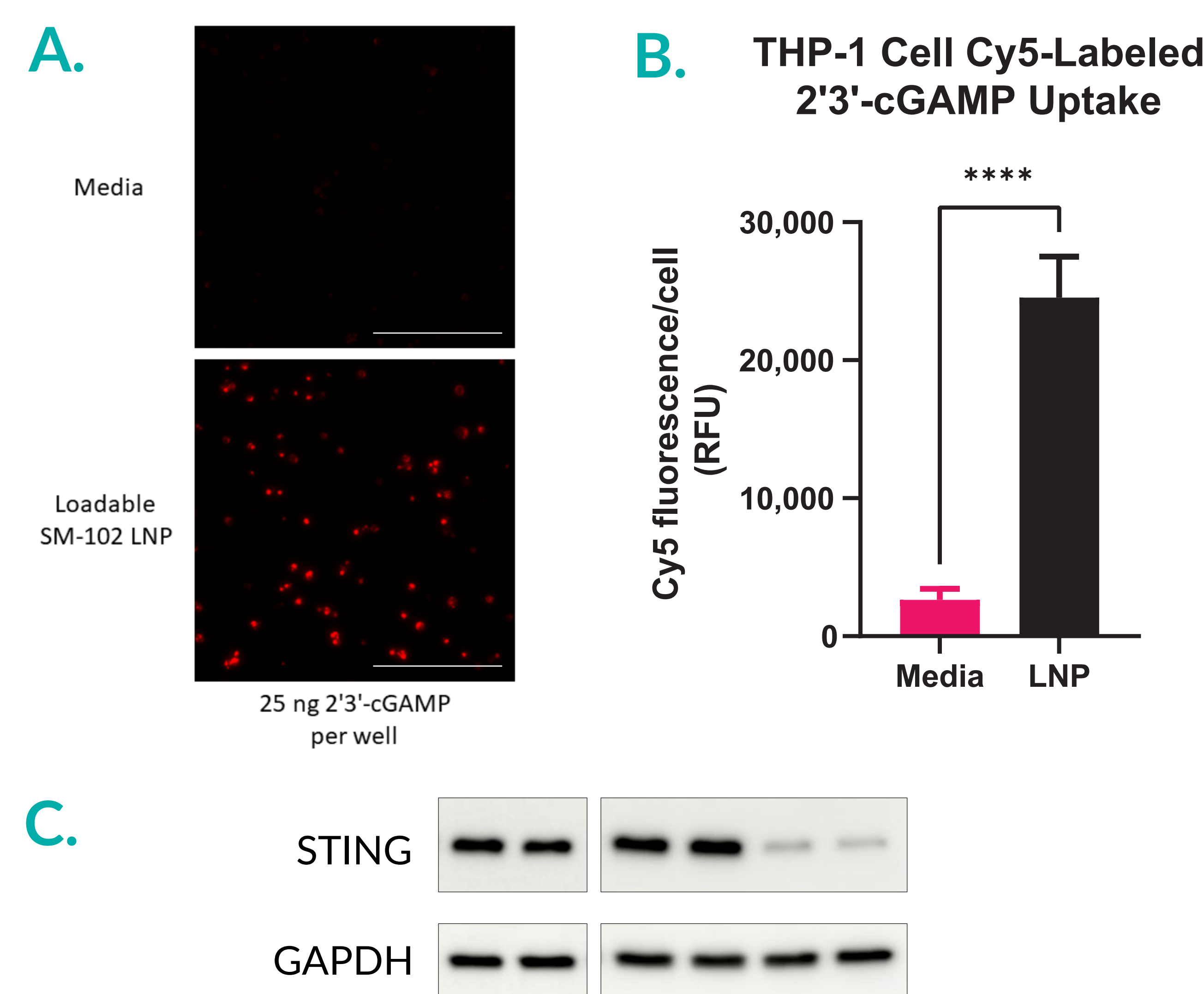


FIGURE 4 – 2'3'-cGAMP levels and cytokine transcript expression in PA-1 cells following siRNA-mediated knockdown of ENPP1.

2'3'-cGAMP released into media from scrambled siRNA-transfected vs. ENPP1 siRNA-transfected PA-1 cells was quantified (A): N=5, ***p<0.001. ENPP1 knockdown led to increased IFNB expression (B) determined via RT-PCR; N=4, *p<0.05.

Preformed loadable SM-102 LNPs enable enhanced 2'3'-cGAMP uptake into THP-1 cells, leading to STING activation



2'3'-cGAMP (μM):	0	0	10	10	10	10
Loadable LNP:	+	+	-	-	+	+

FIGURE 5 – Delivery of 2'3'-cGAMP to THP-1 cells via loadable LNPs leads to STING activation. Uptake of Cy5-2'3'-cGAMP is enhanced when loaded into SM-102 LNPs (A). The quantification is shown in (B). STING is activated in THP-1 cells following treatment with 2'3'-cGAMP loaded LNPs at 10 μM (C). *p<0.05. Bar size = 200 μm.



Download this poster & learn more about our assay kits

RESULTS CONTINUED

PA-1 cells secrete EVs containing ENPP1, which suppressed IL-6 release in 2'3'-cGAMP-stimulated THP-1 macrophages.

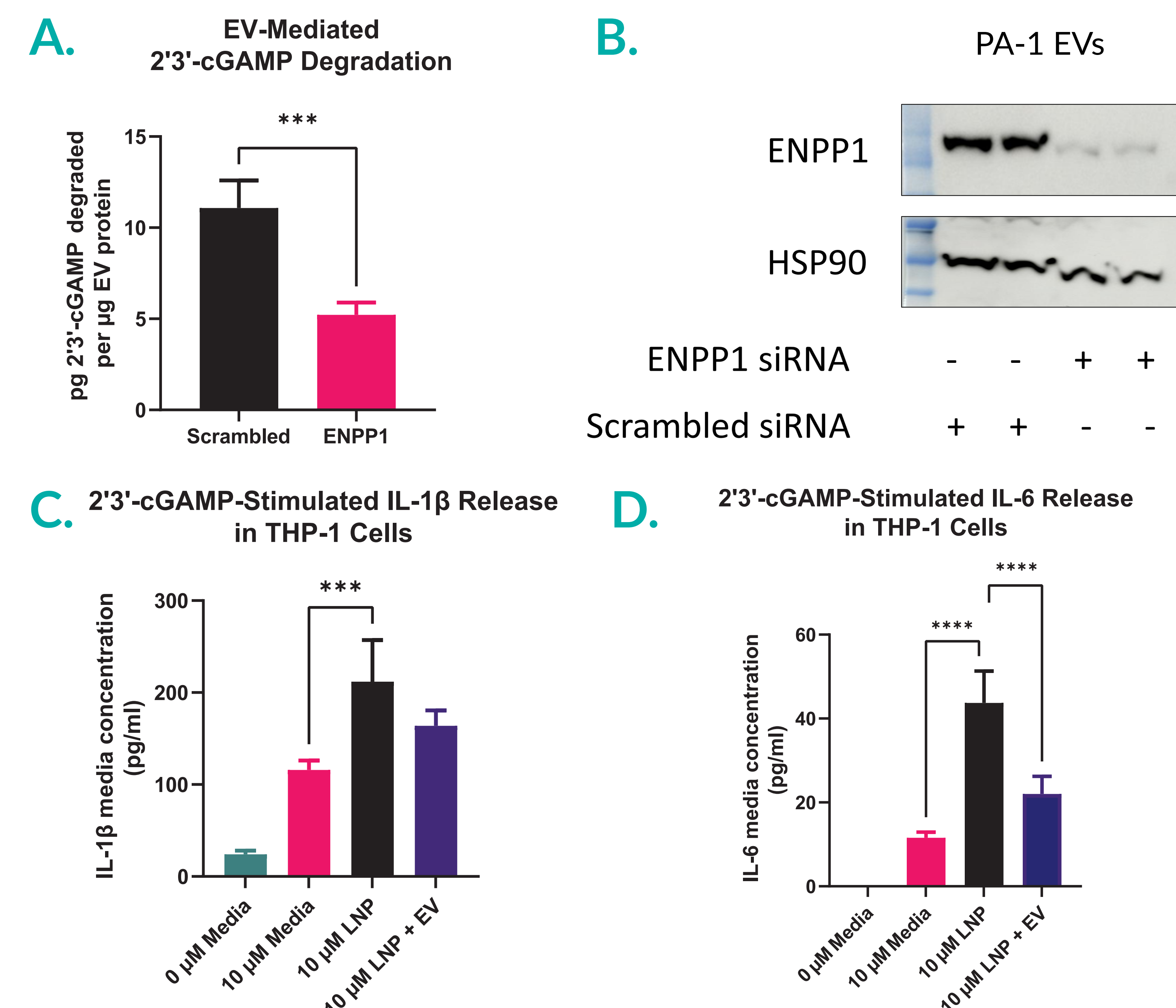


FIGURE 6 – PA-1 cells secrete EVs containing ENPP1, which can modify immune responses.

Degradation of extracellular 2'3'-cGAMP by EVs isolated from scrambled and ENPP1 siRNA-transfected PA-1 cells (A): N=4, ***p<0.001. EV ENPP1 protein levels were reduced following ENPP1 knockdown in PA-1 cells (B). 2'3'-cGAMP stimulated IL-1β and IL-6 production by THP-1 cells was enhanced by LNP-mediated cGAMP delivery (Panels C & D); IL-6 release was blunted by PA-1 cell-derived EVs, N=4, ****p<0.0001.

CONCLUSIONS

- ENPP1 activity was detected in PA-1, Caov-3, and SKOV3 cell lines. Preformed loadable SM-102 LNPs facilitated effective siRNA-mediated ENPP1 knockdown in all three cell lines.
- Among the tested lines, PA-1 cells exhibited the highest ENPP1 activity, which following siRNA-mediated ENPP1 knockdown, led to an increase in extracellular 2'3'-cGAMP levels.
- cGAMP delivery to THP-1 macrophages and activation of STING was enhanced by the use of preformed loadable SM-102 LNPs.
- PA-1 cells secreted EVs containing ENPP1, which blunted 2'3'-cGAMP-stimulated cytokine release in THP-1 macrophages.

References

- Morand, S., Devanaboyina, M., Staats, H., et al. Ovarian cancer immunotherapy and personalized medicine. *Int. J. Mol. Sci.* **22**(12), 6532 (2021).
- Carozza, J.A., Cordova, A.F., Brown, J.A., et al. ENPP1's regulation of extracellular cGAMP is a ubiquitous mechanism of attenuating STING signaling. *Proc. Natl. Acad. Sci. USA* **119**(21), e2119189119 (2022).
- Carozza, J.A., Böhner, V., Nguyen K.C., et al. Extracellular cGAMP is a cancer-cell-produced immunotransmitter involved in radiation-induced anticancer immunity. *Nat. Cancer* **1**(2), 184-196 (2020).
- Li, J., Duran, M.A., Dhanota, N., et al. Metastasis and immune evasion from extracellular cGAMP hydrolysis. *Cancer Discov.* **11**(5), 1212-1227 (2021).
- Ruiz-Fernández de Córdoba, B., Martínez-Monge, R., and Lecanda, F. ENPP1 immunobiology as a therapeutic target. *Clin. Cancer Res.* **29**(12), 2184-2193 (2023).
- Kawaguchi, M., Han, X., Hisada, T., et al. Development of an ENPP1 fluorescence probe for inhibitor screening, cellular imaging, and prognostic assessment of malignant breast cancer. *J. Med. Chem.* **62**(20), 9254-9269 (2019).
- Gangar, M., Goyal, S., Raykar, D., et al. Design, synthesis and biological evaluation studies of novel small molecule ENPP1 inhibitors for cancer immunotherapy. *Bioorg. Chem.* **119**, 105549 (2022).
- Wang, X., Lu, X., Yan, D., et al. Development of novel ecto-nucleotide pyrophosphatase/phosphodiesterase 1 (ENPP1) inhibitors for tumor immunotherapy. *Int. J. Mol. Sci.* **23**(13), 7104 (2022).