



KEY FINDING Differences are observed in the lipid profiles of plasma extracellular vesicles from normal and obese mice.

- EV isolation by precipitation yields a mixture of EV and lipoprotein particles. Efforts are ongoing to improve selectivity of the EV purification.
- DIO mice were found to contain significantly lower protein and lipid levels compared to normal mice.
- Some sphingolipids show the largest differences between healthy and DIO mice. EV lipid profiling reveals significant differences that are not observed in their source plasma.

