

Sponsored by:



Life Sciences

Going Viral

Custom publishing from:

TheScientist

EXPLORING LIFE. INSPIRING INNOVATION.

O.K. Capsid: Where Busy Scientists go to Meet the Virus of Their Dreams

Choosing a viral delivery system is a bit like choosing a partner. Do you share goals? Are you truly compatible? Have you seriously considered asking them to carry a bit of your genetic material? But like any relationship, you need to do your homework before inviting a viral delivery system back to your incubator.

PREFERRED VIRUS #1 ♥♥♥♥♥

ADENOVIRUS
I like big transgenes and I cannot lie.

FRIENDS: 2

- GENE
- HELPER-DEPENDENT ADENOVIRUS

MESSAGES:

HELPER-DEPENDENT ADENOVIRUS: 😞

What's wrong, HDA?

Someone called me gutless today. I don't know why! I'm pretty brave.

Aw, they were just talking about your ability to carry up to 32 kb of foreign DNA!

Enter your message here, please **SEND**

PREFERRED VIRUS #2 ♥♥♥

LENTIVIRUS
I'm a V.I.P. – a very important particle, so I get through the door at all of the best spots.

FRIENDS: 4

- GENE
- FOAMY VIRUS
- RETROVIRUS
- POXVIRUS

MESSAGES:

FOAMY VIRUS: Hey Len, what's your secret to such highly efficient delivery to neurons?

It's simple, really. I just slip in at the nuclear pore!

Enter your message here, please **SEND**

PREFERRED VIRUS #3 ♥♥

ADENO-ASSOCIATED VIRUS
Safety first! Packaging second... or third.

FRIENDS: 4

- GENE
- ADENOVIRUS
- HERPES SIMPLEX VIRUS
- EPSTEIN-BARR VIRUS

MESSAGES:

EPSTEIN-BARR VIRUS: 😞

I'm so tired of all this talk about the hazards of viral gene delivery.

Me too! That's why I'm starting a campaign to limit genome integrations to chromosome 19 where they have no discernable effect!

Enter your message here, please **SEND**

PREFERRED VIRUS #4 ♥

RETROVIRUS
Division is a decision, and I choose dividing cells!

FRIENDS: 4

- GENE
- LENTIVIRUS
- FOAMY VIRUS
- REVERSE TRANSCRIPTASE

MESSAGES:

LENTIVIRUS: You're such a show off! Reverse transcription is all smoke and mirrors.

Oh yeah? Why don't you say hello to my little friend? @RT

REVERSE TRANSCRIPTASE: Hello 🙋

Enter your message here, please **SEND**

Beckman Coulter Life Sciences

Beckman Coulter Life Sciences is dedicated to empowering discovery and scientific breakthroughs. The company's global leadership and world-class service and support deliver sophisticated instrument systems, reagents and services to life science researchers in academic and commercial laboratories, enabling new discoveries in biology-based research and development. A leader in centrifugation and flow cytometry, Beckman Coulter has long been an innovator in particle characterization and laboratory automation, and its products are used at the forefront of important areas of investigation, including genomics and proteomics. For more information, visit beckman.com/home. Follow Beckman Coulter Life Sciences on Twitter @BCLifeSciences; Facebook: BCLifeSciences; and LinkedIn

CENT-2190PST12.16 © 2016 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks used herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries.

Janet looked for a partner that would quickly show her a band.



Her last relationship was plagued by separation anxiety and a lack of clear boundaries. Then Janet met her match. Now she trusts the Optima XPN to meet all her preparative ultracentrifugation needs.

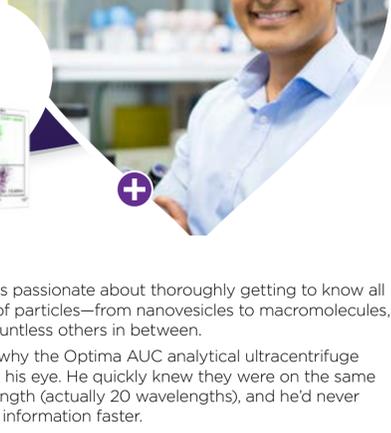
With options ranging from 80,000 – 100,000 RPM, she's not concerned about things going too slowly.

And the Optima XPN's tracking features—as well as multi-level BioSafety* attributes—give her the peace-of-mind she looks for in a long-term relationship.



The Optima XPN ultracentrifuge is ready for any high-performance processing application.

Jay, however, wanted to characterize his options.



Jay was passionate about thoroughly getting to know all types of particles—from nanovesicles to macromolecules, and countless others in between. That's why the Optima AUC analytical ultracentrifuge caught his eye. He quickly knew they were on the same wavelength (actually 20 wavelengths), and he'd never gotten information faster.

But another characterization technique—the CytoFLEX Flow Cytometer**—presented optimal excitation too. Never before had Jay seen a research platform that pushed the boundaries of what's possible with benchtop flow cytometry.

Jay wanted to see as much as possible.

Meet your centrifuge match
Visit info.beckmancoulter.com/thecentrifugeforme to learn more.



Data provided by the Optima AUC (left) and CytoFLEX Flow Cytometer (right).

Going Viral Where Busy Scientists go to Meet the Virus of Their Dreams

