

Show us how much you love your Tecan and win a party for your team!

**WE  OUR
TECAN**



Tecan is proud to have been a true pioneer in the liquid handling market for three decades, and we really hope you value your Tecan as much as we love making them for you.

.....

Send us a photo that captures how important your system is to your team to be in with a chance of winning a huge party for you and 20 of your colleagues worth \$5,000.

The more creative you are the better, as the winning photo will be the one that convinces us your Tecan is the most treasured in the world.

To enter, please go to www.tecan.com/weloveourtecan