



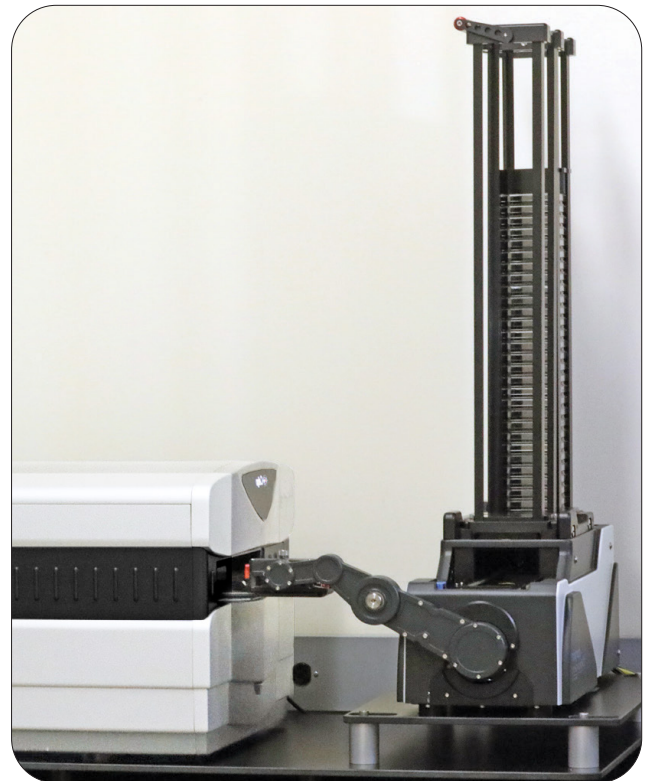
revvity

Take your analysis to the next level: Automation.

Celigo image cytometer plate stacker

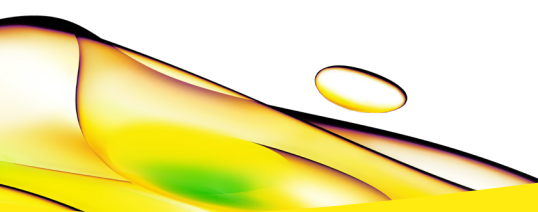
There's no need to "babysit" your plates. With the simple integration of the plate stacker, you are free to work on other tasks while the Celigo™ system screens your stack of plates.

- Up to 50 plate capacity
- 15 sec transfer rate between plates
- Data exports automatically for each plate
- Handles plates with or without lids
- Accommodates from 6-well up to 1536-well plate formats
- Approximate read time for a 96-well plate is under 3.5 minutes.
- Fits on standard lab bench
- Ideal for endpoint assays
- Easy to add on to existing Celigo instruments



| Celigo image cytometer plate stacker

Parameter	Specification
Celigo only dimensions	19 inch wide x 29 inch long x 17 inch high (with plate holder extended)
Celigo + stacker dimensions	25 inch wide x 44 inch long x 40 inch high (Total)
Electrical power	5 Electrical outlets needed at 100-240VAC 50/60Hz (30Amps Total)
Compressor (sold with stacker)	1.2 CFM @ 90 PSI with 1 gallon tank
Operating temperature	15 °C to 25 °C
Operating humidity	10% to 90% RH, non-condensing
Instrument shipping and storage temperature	-18 °C to 65 °C
Instrument shipping and storage humidity	10% to 90% RH, non-condensing



revvity