



How healthy are your cells?

Advanced cell fitness panel for Cellaca[®] MX and Cellometer[®] Spectrum systems.

Multiple assay parameters are used to investigate cell sample fitness. For the first time, these individual tests have been combined into a single kit to provide critical insight into sample health and quality.

1. Viability - AO/PI (acridine orange and propidium iodide)

Proven to be a more accurate cell viability assay compared to trypan blue

2. Vitality/enzymatic activity - Calcein AM/PI

Easily determine whether cells are metabolically active or not

3. Reactive oxygen species - Total RO

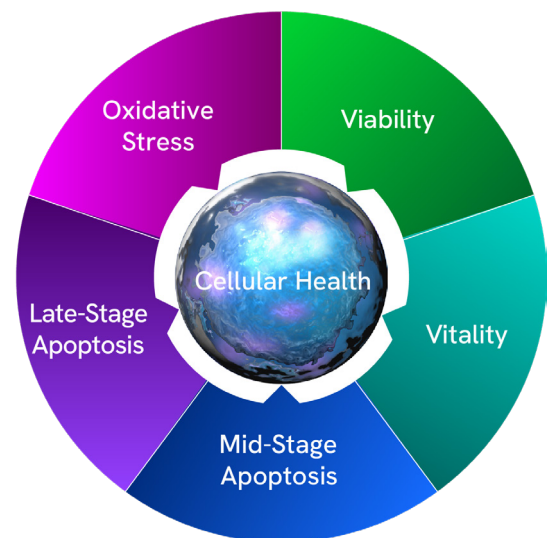
Measure total oxidative stress within your cell culture

4. Mid-stage apoptosis - Annexin V/PI

Rapidly perform a traditional apoptosis assay

5. Late-stage apoptosis - Caspase 3/7

Detect late-stage executioner caspases efficiently



Description	Catalog number	Unit size	Excitation/emission
ViaStain™ cell fitness panel for Cellometer Spectrum	CSK-V0023-1	25 tests per assay	527/655 nm 470/534 nm
ViaStain™ cell fitness panel for Cellaca MX	CSK-V0024-1	1 Cellaca plate per assay	534/655 nm 470/527 nm

Not your average cell counters

Perform cell-based assays and cell counts with one instrument. Both single-sample and high-throughput options are available.

Cellaca MX high-throughput cell counter

High-throughput cell counting

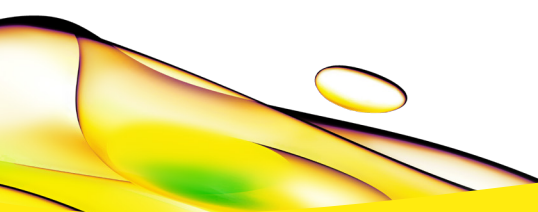
Count 24 plate-based samples in 3 minutes or less



Cellometer Spectrum image cytometry system

Customizable cell counting and image cytometry

Effortlessly run simple two-color assays in 60 seconds



revvity