



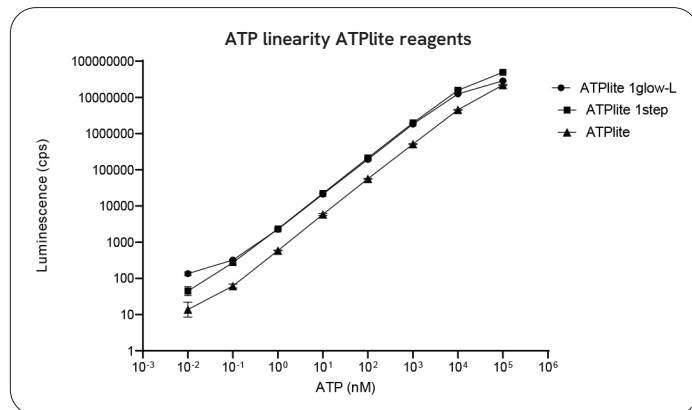
ATPlite - Fast, reliable cytotoxicity and cell proliferation measurement.

Discover a new way forward in measuring cytotoxicity and cell proliferation with our ATPlite™ assay portfolio.

Our ATP luminescence assays offer a robust alternative to traditional colorimetric, fluorometric, and radioisotopic methods for measuring cell viability and proliferation in mammalian cells.

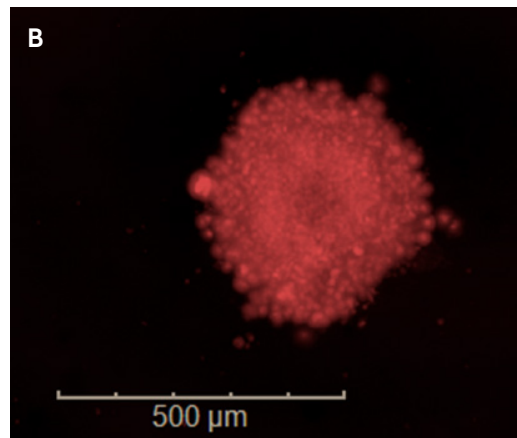
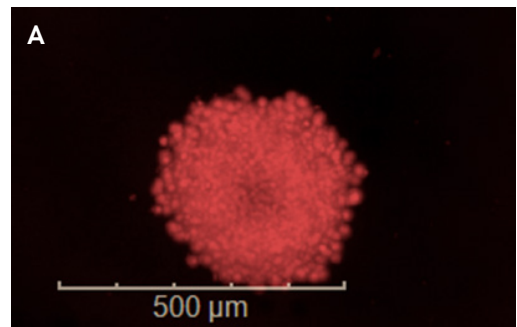
These assays harness firefly luciferase to detect ATP, producing light in direct proportion to ATP concentration—a reliable marker of metabolically active cells.

Explore our full range of 2D and 3D cytotoxicity and cell proliferation assays to find the ideal solution for your research need.



ATPlite kits standard curves.

DRAQ7 staining 3D Assay spheroid permeabilization



ATPlite 1glow-F 3D assay

LMSU gastric adherent cell line spheroids grown 7 days and treated with ATPlite 1glow-F (A) and a competitive reagent (B). DRAQ7 staining measures permeabilization and shows ATPlite 1glow-F penetrates the spheroid.

Choose the right format for your needs:

- **ATPlite 1glow** - Single addition with extended signal stability (> 3 hours), wide dynamic range, and high-throughput automation friendly. Available as ATPlite 1glow-F, frozen ready-to-use and ATPlite 1glow-L lyophilized formats for 2D and 3D cell cultures.
- **ATPlite 1step** - Single addition assay for streamlined workflows
- **ATPlite** - Extended signal stability for flexible timing
- **3D applications** - Both ATPlite 1step and ATPlite are available in specialized formats for 3D spheroids, microtissues, and organoids—providing more physiologically relevant results that better bridge the gap between traditional 2D cultures and in vivo models.

ATPlite cytotoxicity and cell proliferation assays: choose the right detection reagents based on your application

Product	Signal half-life	Kit size	No. of data points			Part No.
			96-w	384-w	1536-w	
ATPlite 1step	30m	10 mL	100	400	3,300	6016736
		100 mL	1,000	4,000	33,000	6016731
		1,000 mL	10,000	40,000	333,000	6016739
ATPlite	5h	300 assays	300	1,200	9,900	6016943
		1,000 assays	1,000	4,000	33,000	6016941
		5,000 assays	5,000	20,000	165,000	6016947
		10,000 assays	10,000	40,000	333,000	6016949
ATPlite 1glow-L	> 3h	10 mL	100	400	3,300	6096746
		100 mL	1,000	4,000	33,000	6096741
		1000 mL	10,000	40,000	333,000	6096749
ATPlite 1glow-F	> 3h	10 mL	100	400	3,300	6096736
		100 mL	1,000	4,000	33,000	6096731
		450 mL	4,500	18,000	112,000	6096739
ATPlite 1step 3D	30m	100 assays	100	400	3,300	6066736
		1000 assays	1,000	4,000	33,000	6066731
ATPlite 3D	5h	300 assays	300	1,200	9,900	6066943

- All lyophilized kits include at least one ATP standard vial
- ATPlite 1step 10 mL includes a 96-well CulturPlate™ with a lid
- The 10mL ATPlite 1step 3D and ATPlite 3D kits include a CellCarrier™ Spheroid ULA 96-well microplate, an OptiPlate™, 96-well, white microplate, and 4 TopSeal™-A PLUS clear adhesive microplate seals

Complementary Revvity products	Part no.
ATP Standard	6016841
CellCarrier Spheroid ULA 96-well microplates	6055330 (10 lidded, individually-wrapped plates)
	6055334 (40 lidded plates - two bags with 20/bag)
OptiPlate™, 96-well, white	6055290 (case of 50)
	6055299 (case of 200)
OptiPlate, 384-well, white	6007290 (case of 50)
	6007299 (case of 200)
AlphaPlate™ 384-well, gray	6057350 (case of 50)
	6057359 (case of 200)
TopSeal™-A PLUS microplate seals	6050185 (box of 100)



revvity