



CELL HEALTH ASSAYS

Product	What is Being Measured?	Cells Detected			Signal Detected				Compatible Plate Format			Minimum Incubation Time	96-Well Sensitivity	Cat.#
		Live Cells	Dead Cells	Apoptotic Cells	Luminescence	Fluorescence	Color Absorbance	Real-Time Detection	96	384	1536			
RealTime-Glo™ MT Cell Viability Assay	Active metabolism	✓			✓			✓	✓	✓	✓	Real-time monitoring	<100 cells/well	G9711
CellTiter-Glo® 2.0 Assay	ATP from viable cells	✓			✓				✓	✓	✓	10 minutes	15 living cells	G9241
CellTiter-Glo® Luminescent Cell Viability Assay		✓			✓				✓	✓	✓	10 minutes	10 living cells	G7570
CellTiter-Glo® 3D Cell Viability Assay		✓			✓				✓	✓	✓	30 minutes		G9681
CellTiter-Glo® One Solution Assay		✓			✓				✓	✓	✓	10 minutes		G8461
CellTiter-Fluor™ Cell Viability Assay		Retained live-cell protease activity	✓				AFC 400 _{Ex} /505 _{Em}			✓	✓	✓	30 minutes	40 living cells
CellTiter-Blue® Cell Viability Assay	Resazurin reduction	✓				Resofurin 560 _{Ex} /590 _{Em}	570nm		✓	✓		1–4 hours	400 living cells	G8080
CellTiter 96® Non-Radioactive Cell Proliferation Assay	MTT reduction	✓					570nm		✓			4 hours		G4000
CellTiter 96® Aqueous One Solution Cell Proliferation Assay	MTS reduction	✓					490nm		✓	✓		1–4 hours	1,000 living cells	G3582
CellTiter 96® Aqueous MTS Reagent Powder		✓					490nm		✓	✓		1–4 hours	1,000 living cells	G1112
MultiTox-Fluor Multiplex Cytotoxicity Assay	Live- and dead-cell protease activity	✓	✓			AFC 400 _{Ex} /505 _{Em} R110 485 _{Ex} /520 _{Em}			✓	✓	✓	30 minutes	40 living cells, 10 dead cells	G9200
MultiTox-Glo Multiplex Cytotoxicity Assay		✓	✓		✓	AFC 400 _{Ex} /505 _{Em}			✓	✓	✓	30+15 minutes	40 living cells, 10 dead cells	G9270
LDH-Glo™ Cytotoxicity Assay	LDH release		✓		✓				✓	✓	✓	30 minutes– 1 hour	10 dead cells	J2380
CytoTox-ONE™ Homogeneous Membrane Integrity Assay			✓				Resofurin 560 _{Ex} /590 _{Em}			✓	✓	10 minutes	200 dead cells	G7890
CytoTox 96® Non-Radioactive Cytotoxicity Assay				✓				490nm		✓	✓	30 minutes		G1780
CytoTox-Fluor™ Cytotoxicity Assay		Released dead-cell protease activity		✓			R110 485 _{Ex} /520 _{Em}			✓	✓		30 minutes	10 dead cells
CytoTox-Glo™ Cytotoxicity Assay				✓		✓				✓	✓	✓	15 minutes	10 dead cells
CellTox™ Green Cytotoxicity Assay	DNA staining upon membrane disruption		✓			480-500 _{Ex} / 520-530 _{Em}		✓	✓	✓	✓	15 minutes		G8741
ApoTox-Glo™ Triplex Assay	Live-cell protease + Dead-cell protease + Caspase 3/7 activity	✓	✓	✓	✓	AFC 400 _{Ex} /505 _{Em} R110 485 _{Ex} /520 _{Em}			✓	✓		30+30 minutes	40 living cells, 100 apoptotic cells	G6320
RealTime-Glo™ Annexin V Apoptosis and Necrosis Assay	Annexin HiBIT luciferase binding to phosphatidylserine + DNA staining upon membrane disruption		✓	✓	✓	✓		✓	✓	✓	✓	Real-time monitoring	100 apoptotic cells	JA1011
ApoLive-Glo™ Multiplex Assay	Live-cell protease + Caspase 3/7 activity	✓		✓	✓	AFC 400 _{Ex} /505 _{Em}			✓	✓		30+30 minutes	40 living cells, 100 apoptotic cells	G6410
Apo-ONE® Homogeneous Caspase-3/7 Assay	Caspase-3/7 activity			✓		R110 499 _{Ex} /521 _{Em}			✓	✓	✓	30 minutes	625 apoptotic cells	G7790
Caspase-Glo® 3/7 Assay					✓	✓				✓	✓	✓	20 minutes	100 apoptotic cells
Caspase-Glo® 8 Assay	Caspase-8 activity			✓	✓				✓	✓		20 minutes	1,000 apoptotic cells	G8200
Caspase-Glo® 9 Assay	Caspase-9 activity			✓	✓				✓	✓		20 minutes	1,500 apoptotic cells	G8210