



Assays for **GLOMAX®-96 MICROPLATE LUMINOMETER**



DUAL-REPORTER ASSAYS

Dual-Glo™ Luciferase Assay	Dual Injector System E2920
Dual-Luciferase® Reporter Assay System	Dual Injector System E1910

FIREFLY LUCIFERASE REPORTER ASSAYS

Luciferase Assay System	E1500
Bright-Glo™ Luciferase Assay System	E2610
Steady-Glo® Luciferase Assay System	E2510
ONE-Glo® Luciferase Assay System	E6110

RENILLA LUCIFERASE REPORTER ASSAYS

EnduRen™ Live Cell Substrate	E6481
ViviRen™ Live Cell Substrate	E6491
Renilla Luciferase Assay System	E2810
Renilla Glo	E2720

CELL VIABILITY ASSAYS

CellTiter-Glo® Luminescent Cell Viability Assay	G7570
GSH/GSSG-Glo™ Assay	V6611
GSH-Glo™ Glutathione Assay	V6911
BacTiter-Glo™ Microbial Cell Viability Assay	G8320
CytoTox-Glo™ Cytotoxicity Assay	G9290

APOPTOSIS ASSAYS

Caspase-Glo® 3/7 Assay	G8091
Caspase-Glo® 8 Assay	G8201
Caspase-Glo® 9 Assay	G8211
Caspase-Glo® 2 Assay	G0940
Caspase-Glo® Assay	G0970
cAMP-GLO™ Max Assay	V1681

CELL SIGNALING

cAMP-Glo™ Assay	V1501
PDE-Glo™ Phosphodiesterase Assay	V1361

CYP450 SCREENING SYSTEMS

P450-Glo™ CYP1A2 Screening System	V9770
P450-Glo™ CYP2C9 Screening System	V9790
P450-Glo™ CYP3A4 Screening System	V9800
P450-Glo™ CYP2C19 Screening System	V9880
P450-Glo™ CYP2D6 Screening System	V9890
P450-Glo™ CYP3A4 Screening System (Luciferin - PPXE) DMSO-Tolerant Assay	V9910



Assays for
**GLOMAX®-96 MICROPLATE
LUMINOMETER**

LUMINESCENT ENZYME ASSAYS

UGT-Glo™ Assay	V2081
MAO-Glo™ Assay	V1401
Pgp-Glo™ Assay	V3591

LUMINESCENT KINASE ASSAYS

ADP-Glo™ Kinase Assay	V9101
Kinase-Glo® Luminescent Kinase Assay	V6711
Kinase-Glo® Plus Luminescent Kinase Assay	V3771
Kinase-Glo® Max Luminescent Kinase Assay	V6071

PROTEASE ASSAYS

DPPIV-Glo™ Protease Assay	G8350
Calpain-Glo™ Protease Assay	G8501
Protease-Glo Assay	G9451
Proteasome-Glo™ Chymotrypsin-Like Assay	G8621
Proteasome-Glo™ Trypsin-Like Assay	G8631
Proteasome-Glo™ Caspase-Like Assay	G8641
Proteasome-Glo™ 3-Substrate System	G8531
Dub-Glo™ Protease Assay	G6260
Protease-Glo™ Assay	G9451

