

<b>Cat #</b>	<b>Cell Line</b>
EL-001	Ba/F3 Control Cell Line EL-001
EL-002	EL-002 EGFR WT Ba/F3 Stable Cell Line
EL-003	EL-003 EGFR (L858R) Ba/F3 Stable Cell Line
EL-004	EL-004 EGFR (L858R+T790M) Ba/F3 Stable Cell Line
EL-005	EL-005 EGFR (Del19 (Del746-750)) HCC827 Stable Cell Line
EL-006	EL-006 EGFR (Del19 (Del746-750)+T790M) HCC827 Stable Cell Line
EL-007	Control (Empty Vector) HCC827 Stable Cell Line EL-007
EL-008	EL-008 EGFR( D770_N771insSVD ) BaF3 Stable Cell Line
EL-009	EL-009 EGFR (A763_Y764insFQEA) BaF3 Stable Cell Line
EL-010	EL-010 EGFR (A767_dupASV) BaF3 Stable Cell Line
EL-011	EL-011 EGFR (Del19 (del747-753InsS)) BaF3 Stable Cell Line
EL-014	EL-014 EGFR (Del19 (del747-753InsS)-T790M) BaF3 Stable Cell Line
EL-016	EL-016 EGFR (L858R) CHO Stable Cell Line
EL-017	EL-017 EGFR (T790M) CHO Stable Cell Line
EL-018	EL-018 EGFR (M766_A767insAI) BaF3 Stable Cell Line
EL-019	EL-019 EGFR (H773_V774insH) BaF3 Stable Cell Line
EL-020	EL-020 EGFR (Y764_V765insHH) BaF3 Stable Cell Line
EL-021	EL-021 EGFR (Del19 (Del746-750)) BaF3 Stable Cell Line
EL-024	EL-024 EGFR (Del19 (Del746-750) + T790M) BaF3 Stable Cell Line
EL-031	HER2 (WT) BaF3 Stable Cell Line EL-0031
EL-032	HER2 (S310F) BaF3 Stable Cell Line EL-0032
EL-033	HER2 (A775_G776insYVMA) BaF3 Stable Cell Line EL-0033
EL-034	HER2 (P780_Y781insGSP) BaF3 Stable Cell Line EL-0034
EL-035	HER2 (P95) BaF3 Stable Cell Line EL-0035
EL-037	EL-037 EGFR (A767_dupASV, C797S) BaF3 Stable Cell Line
EL-038	EL-038 EGFR (D770_N771insSVD, C797S) BaF3 Stable Cell Line
EL-039	EL-039 EGFR (H773_V774insH, C797S) BaF3 Stable Cell Line
EL-040	EL-040 EGFR (L718Q, T790M, L858R) BaF3 Stable Cell Line
EL-042	EL-042 EGFR (T790M, L792H, L858R) BaF3 Stable Cell Line
EL-043	EL-043 EGFR (L792H, L858R) BaF3 Stable Cell Line
EL-044	HER2 (A775_G776insYVMA, C805S) BaF3 Cells EL-0044
EL-046	EL-046 EGFR (Del19 (Del746-750) + T790M) and cMET BaF3 Stable Cell
EL-1001	GSK3B Hela Stable Cell Line EL-1001
EL-1002	GSK3B HEK293 Stable Cell Line EL-1002
EL-1003	MMP9 HEK293 Stable Cell Line EL-1003
PC-001	Hela Cells
PC-002	HEK293
PC-003	A549 Cells
PC-004	NIH/3T3
PC-005	HepG2

PC-006	Jurkat Cells
PC-007	K562 Cells
PC-008	MCF7
PC-010	COS-7
PC-011	SF9
SL-0001	NFkB Luciferase Reporter HeLa Stable Cell Line SL-0001
SL-0002	Estrogen Receptor Luciferase Reporter T47D Stable Cell Line SL-0002
SL-0003	Stat3 Luciferase Reporter Hela Stable Cell Line SL-0003
SL-0004	Stat1 Luciferase Reporter Hela Stable Cell Line SL-0004
SL-0005	HIF Luciferase Reporter NIH3T3 Stable Cell Line SL-0005
SL-0006	NFkB Luciferase Reporter NIH 3T3 Stable Cell Line SL-0006
SL-0007	p53 Luciferase Reporter RKO Stable Cell Line SL-0007
SL-0008	AR Luciferase Reporter MDA-MB-453 Stable Cell Line SL-0008
SL-0009	GR Luciferase Reporter MDA-MB-453 Stable Cell Line SL-0009
SL-0010	NRF2/ARE Luciferase Reporter MCF7 Stable Cell Line SL-0010
SL-0011	p53 Luciferase Reporter Hela Stable Cell Line SL-0011
SL-0012	NFkB Luciferase Reporter HEK293 Stable Cell Line SL-0012
SL-0013	NFkB Luciferase Reporter MCF7 Stable Cell Line SL-0013
SL-0014	NFkB Luciferase Reporter A549 Stable Cell Line SL-0014
SL-0015	TCF/LEF Luciferase Reporter HEK293 Stable Cell Line SL-0015
SL-0016	SMAD/TGFbeta Luciferase Reporter HepG2 Stable Cell Line SL-0016
SL-0017	NFkB Luciferase Reporter HepG2 Stable Cell Line SL-0017
SL-0018	NFAT Luciferase Reporter Hela Stable Cell Line SL-0018
SL-0019	AP-1 Luciferase Reporter Hela Stable Cell Line SL-0019
SL-0020	CREB Luciferase Reporter HEK293 Stable Cell Line SL-0020
SL-0021	Glucocorticoid Receptor Luciferase Reporter Hela Stable Cell Line SL-0021
SL-0022	TCF/LEF Luciferase Reporter Hela Stable Cell Line SL-0022
SL-0023	HIF Luciferase Reporter Hela Stable Cell Line SL-0023
SL-0024	ATF6 Luciferase Reporter CHO-K1 Stable Cell Line SL-0024
SL-0025	CHOP Luciferase Reporter Mia_Paca2 Stable Cell Line SL-0025
SL-0026	NFkB Luciferase Reporter Neuro2a Stable Cell Line SL-0026
SL-0027	HIF Luciferase Reporter Neuro2a Stable Cell Line SL-0027
SL-0028	TCF/LEF Luciferase Reporter CHO-K1 Stable Cell Line SL-0028
SL-0029	NFAT Luciferase Reporter NIH 3T3 Stable Cell Line SL-0029
SL-0030	SMAD/TGFbeta Luciferase Reporter NIH/3T3 Stable Cell Line SL-0030
SL-0031	CREB Luciferase Reporter NIH/3T3 Stable Cell Line SL-0031
SL-0032	NFAT Luciferase Reporter Jurkat Stable Cell Line SL-0032
SL-0033	NFkB Luciferase Reporter MEF Stable Cell Line SL-0033
SL-0034	HIF Luciferase Reporter Cos-7 Stable Cell Line SL-0034
SL-0035	IRF Luciferase Reporter HEK293 Stable Cell Line SL-0035
SL-0038	Control Hela Luciferase Reporter Stable Cell Line SL-0038

SL-0039	Control HepG2 Luciferase Reporter Stable Cell Line SL-0039
SL-0040	ELK-TAD Luciferase Reporter HEK293 Stable Cell Line SL-0040
SL-0041	ELK-TAD Luciferase Reporter Hela Stable Cell Line SL-0041
SL-0042	NRF2/ARE Luciferase Reporter HEK293 Stable Cell Line SL-0042
SL-0043	NFkB Luciferase Reporter MDA-MB-231 Stable Cell Line SL-0043
SL-0046	NRF2/ARE Luciferase Reporter HepG2 Stable Cell Line SL-0046
SL-0047	NRF2/ARE Luciferase Reporter NIH3T3 Stable Cell Line SL-0047
SL-0048	IL-6 Responsive Luciferase Reporter HEK293 Cell Line SL-0048
SL-0049	IRF Luciferase Reporter HepG2 Stable Cell Line SL-0049
SL-0050	NFkB Luciferase Reporter Jurkat Stable Cell Line SL-0050
SL-0051	SMAD/BMP Responsive Luciferase Reporter HEK293 Stable Cell Line SL-0051
SL-0052	IFNa-Luciferase Reporter HeLa Stable Cell Line SL-0052
SL-0053	MRF Luciferase Reporter HEK293 Stable Cell Line SL-0053
SL-0055	FXR Luciferase Reporter HepG2 Stable Cell Line SL-0055
SL-0065	Amino acid deprivation-AARE Luciferase Reporter MEF Stable Cell Line SL-0065
SL-0066	Amino acid deprivation-AARE Luciferase Reporter 293 Stable Cell Line SL-0066
SL-0067	ER Stress-ERSE Luciferase Reporter MEF Stable Cell Line SL-0067
SL-0068	ER Stress-ERSE Luciferase Reporter HEK293 Stable Cell Line SL-0068
SL-0069	ER Stress-ERSE Luciferase Reporter HepG2 Stable Cell Line SL-0069
SL-0071	Stat3 Luciferase Reporter HEK293 Stable Cell Line SL-0071
SL-0072	p53 Luciferase Reporter HCT116 Stable Cell Line SL-0072
SL-0074	GAS/Stat1 Luciferase Reporter HCT116 Stable Cell Line SL-0074
SL-0075	XRE Luciferase Reporter HEK293 Stable Cell Line SL-0075
SL-0077	ELK Luciferase Reporter HEK293 Stable Cell Line SL-0077
SL-0078	NFAT Luciferase Reporter HEK293 Stable Cell Line SL-0078
SL-0079	ATF4 Luciferase Reporter HepG2 Stable Cell Line SL-0079
SL-0080	ATF4 Luciferase Reporter HEK293 Stable Cell Line SL-0080
SL-0081	Stat3 Luciferase Reporter HepG2 Stable Cell Line SL-0081
SL-0082	ATF6 Luciferase Reporter HepG2 Stable Cell Line SL-0082
SL-0083	CHOP Luciferase Reporter HEK293 Stable Cell Line SL-0083
SL-0084	ATF6 Luciferase Reporter HEK293 Stable Cell Line SL-0084
SL-0085	AP1 Luciferase Reporter HEK293 Stable Cell Line SL-0085
SL-0086	RAR Luciferase Reporter HEK293 Stable Cell Line SL-0086
SL-0100	p53 Stable Cell Line SL-0100
SL-0101	NFkB (p65) Stable Cell Line SL-0101
SL-0102	Luciferase HeLa Stable Cell Line SL-0102
SL-0103	Luciferase HEK293 Stable Cell Line SL-0103
SL-0104	Luciferase A549 Stable Cell Line SL-0104
SL-0113	FLT3 D835Y Ba/F3 Stable Cell Line SL-0113
SL-0114	FLT3 D835V Ba/F3 Stable Cell Line SL-0114
SL-0115	FLT3 D835E Ba/F3 Stable Cell Line SL-0115

SL-0116	FLT3 D835N Ba/F3 Stable Cell Line SL-0116
SL-0117	FLT3 D835F Ba/F3 Stable Cell Line SL-0117
SL-0118	FLT3 D835H Ba/F3 Stable Cell Line SL-0118
SL-0119	FLT3 WT Ba/F3 Stable Cell Line SL-0119
SL-0200	Jak2WT Expressing Ba/F3 Cell Line
SL-0201	Jak2 V617F Expressing Ba/F3 Cell Line
SL-0202	Jak2 V617F + EPOR expressing Ba/F3 cell line SL-0202 (G418 resistance)
SL-3000	UAS Luciferase Reporter HEK293 Stable Cell Line
SL-3001	GR LBD-driven GAL4 reporter HEK 293 Stable Cell Line SL-3001
SL-3001- HS	Ultra Sensitive GR LBD-driven GAL4 reporter HEK293 stable cell line
SL-3001- HS	Ultra Sensitive GR LBD-driven GAL4 reporter HEK293 stable cell line
SL-3002	PPAR-gamma LBD-Driven GAL4 Reporter HEK 293 Stable Cell Line SL-3002
SL-3002- HS	Ultra Sensitive PPAR $\gamma$ LBD-Driven GAL4 Reporter HEK293 Stable Cell Line
SL-3003	PPAR-alpha LBD-Driven GAL4 Reporter HEK 293 Stable Cell Line SL-3003
SL-3003- HS	Ultra Sensitive PPAR $\alpha$ LBD-Driven GAL4 Reporter HEK293 Stable Cell Line
SL-3004	PPAR-delta LBD-Driven GAL4 Reporter HEK 293 Stable Cell Line SL-3004
SL-3004- HS	Ultra Sensitive PPAR $\delta$ LBD-Driven GAL4 Reporter HEK293 Stable Cell Line
SL-3005	Ultra Sensitive PXR LBD-Driven GAL4 Reporter HEK 293 Stable Cell Line
SL-4000	UAS Luciferase Reporter HeLa Stable Cell Line
SL-4001	NFKB Secreted Luciferase Reporter HeLa Stable Cell Line SL-4001
SL-4012	NFKB Secreted Luciferase Reporter HEK293 Stable Cell Line SL-4012
SL-5000	UAS Luciferase Reporter CHO Stable Cell Line