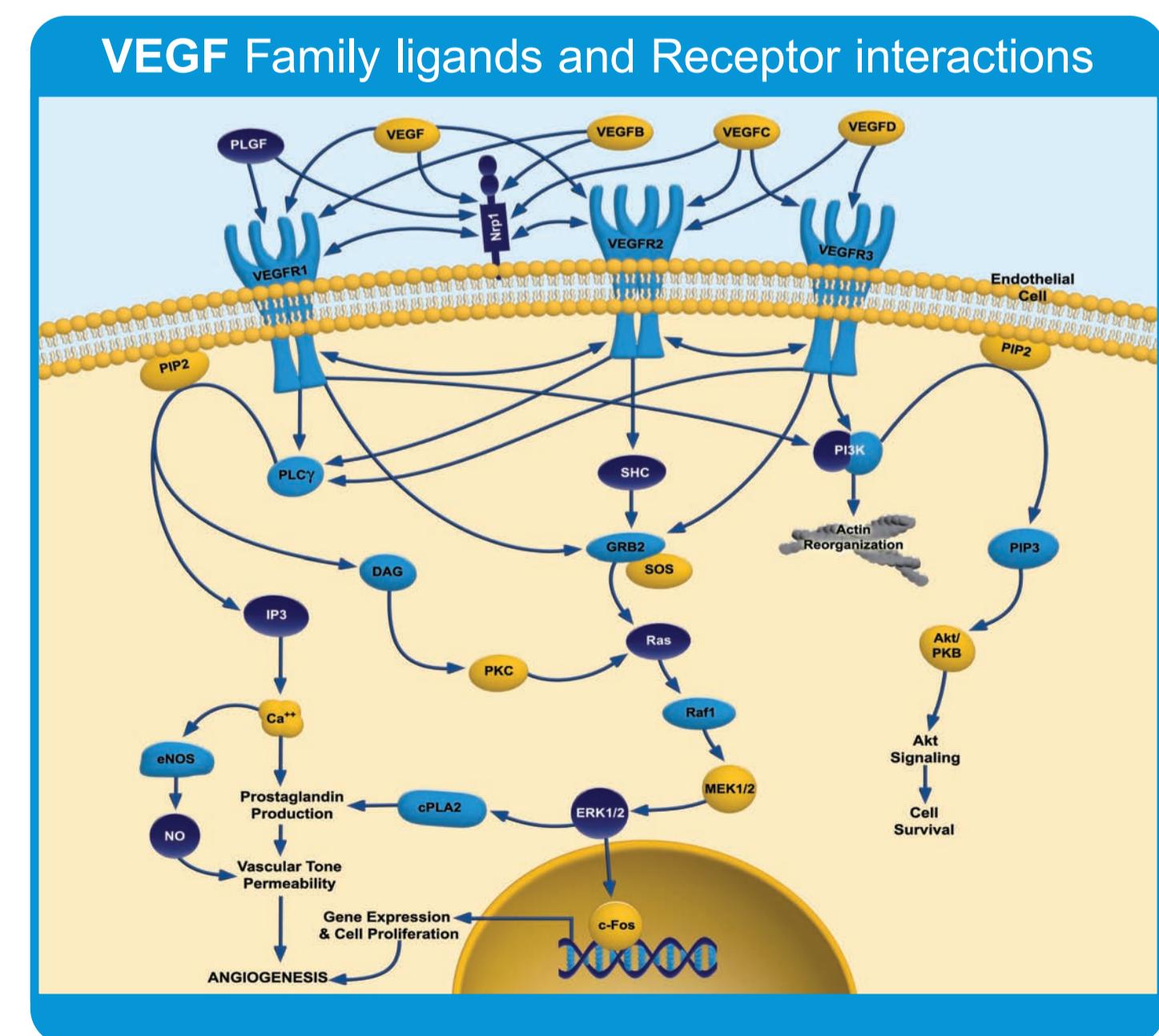
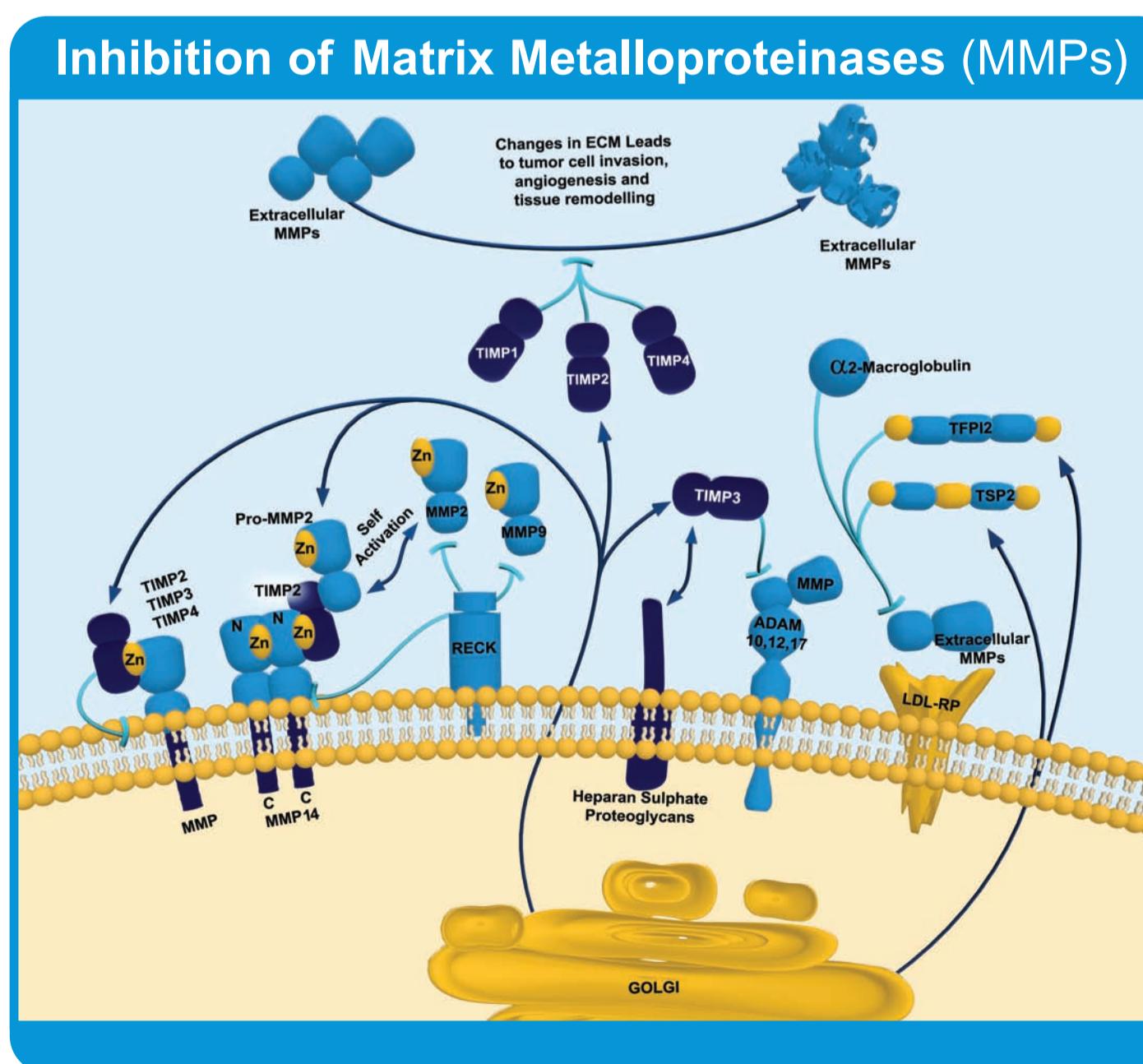
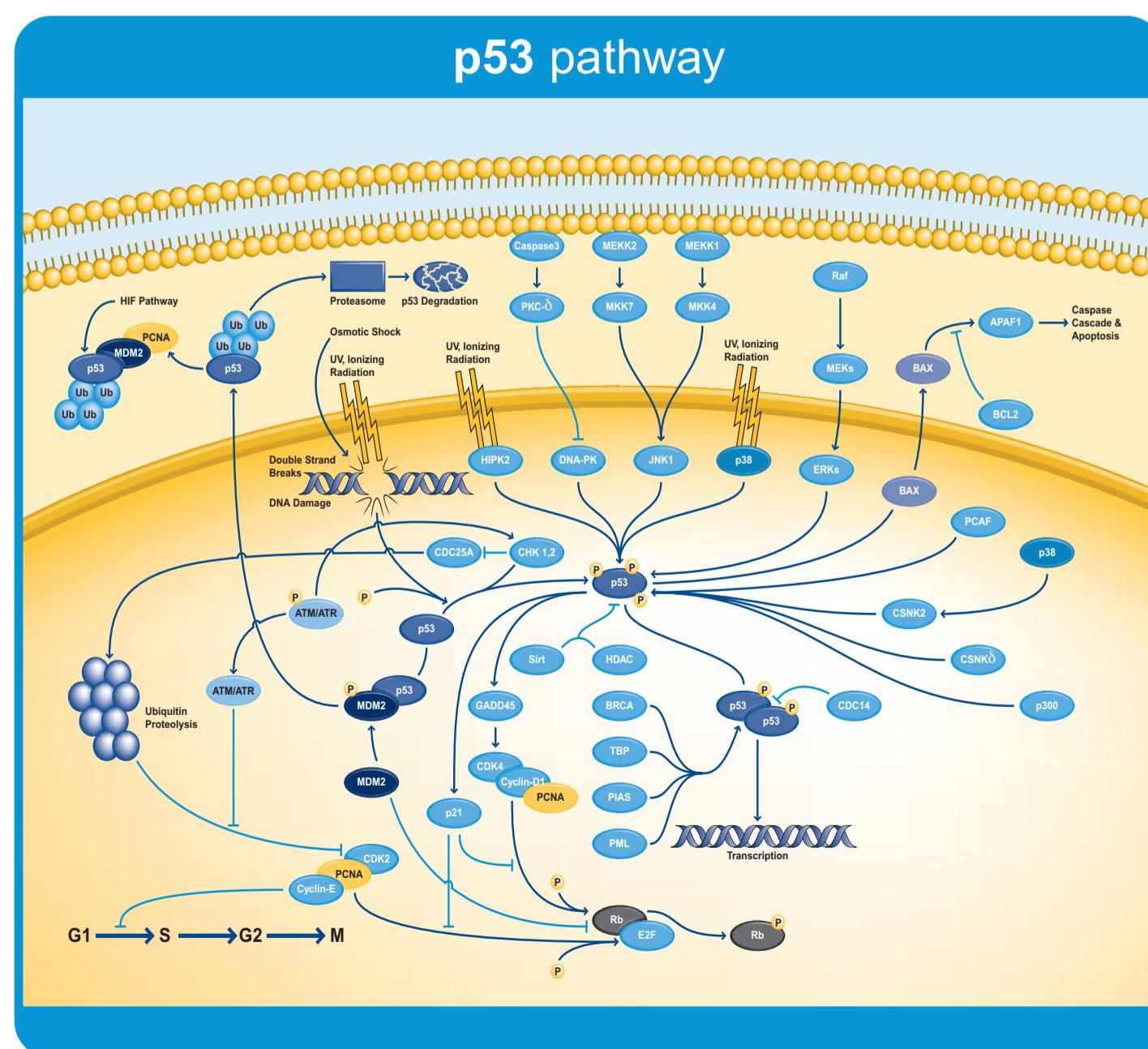
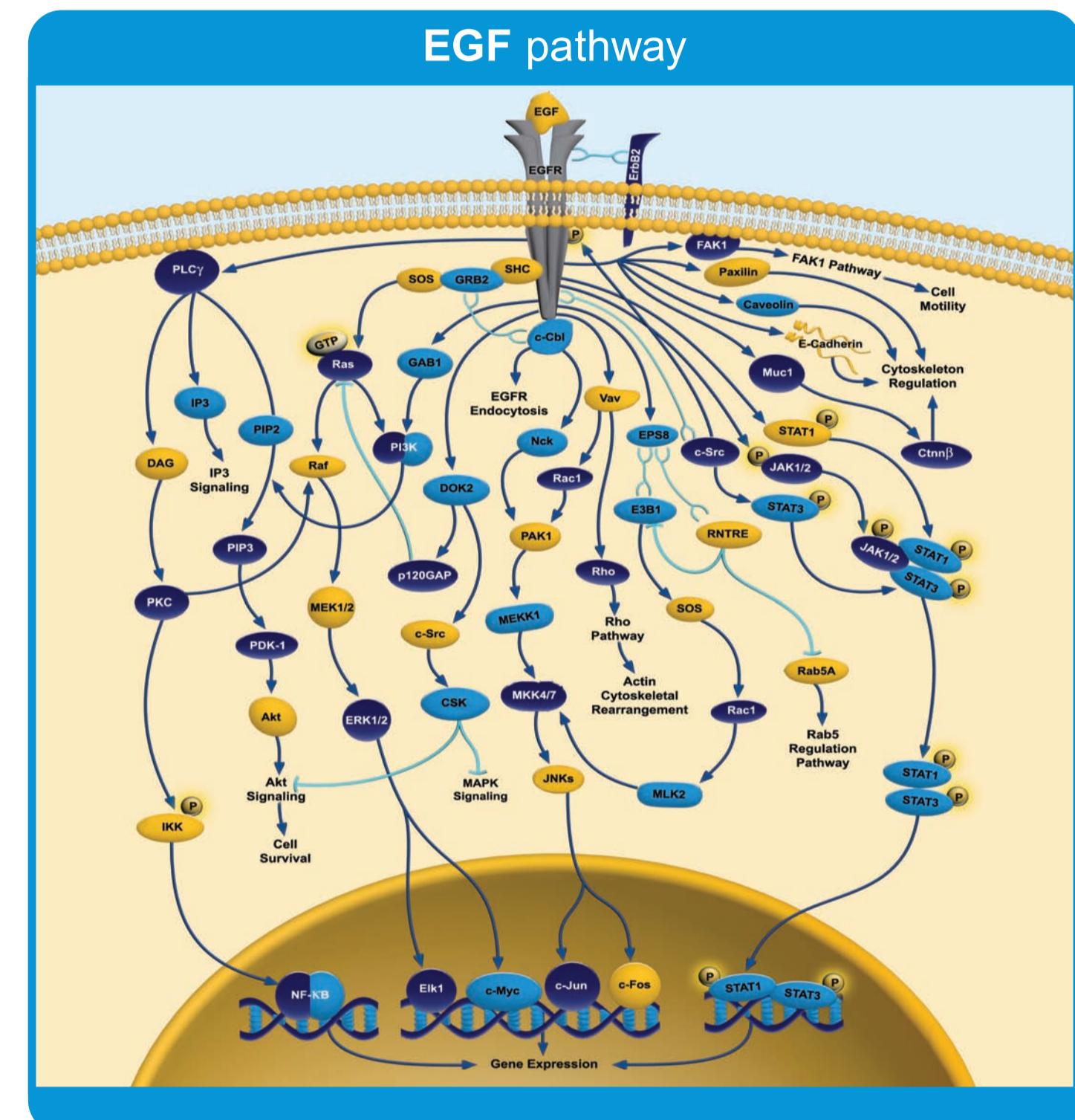
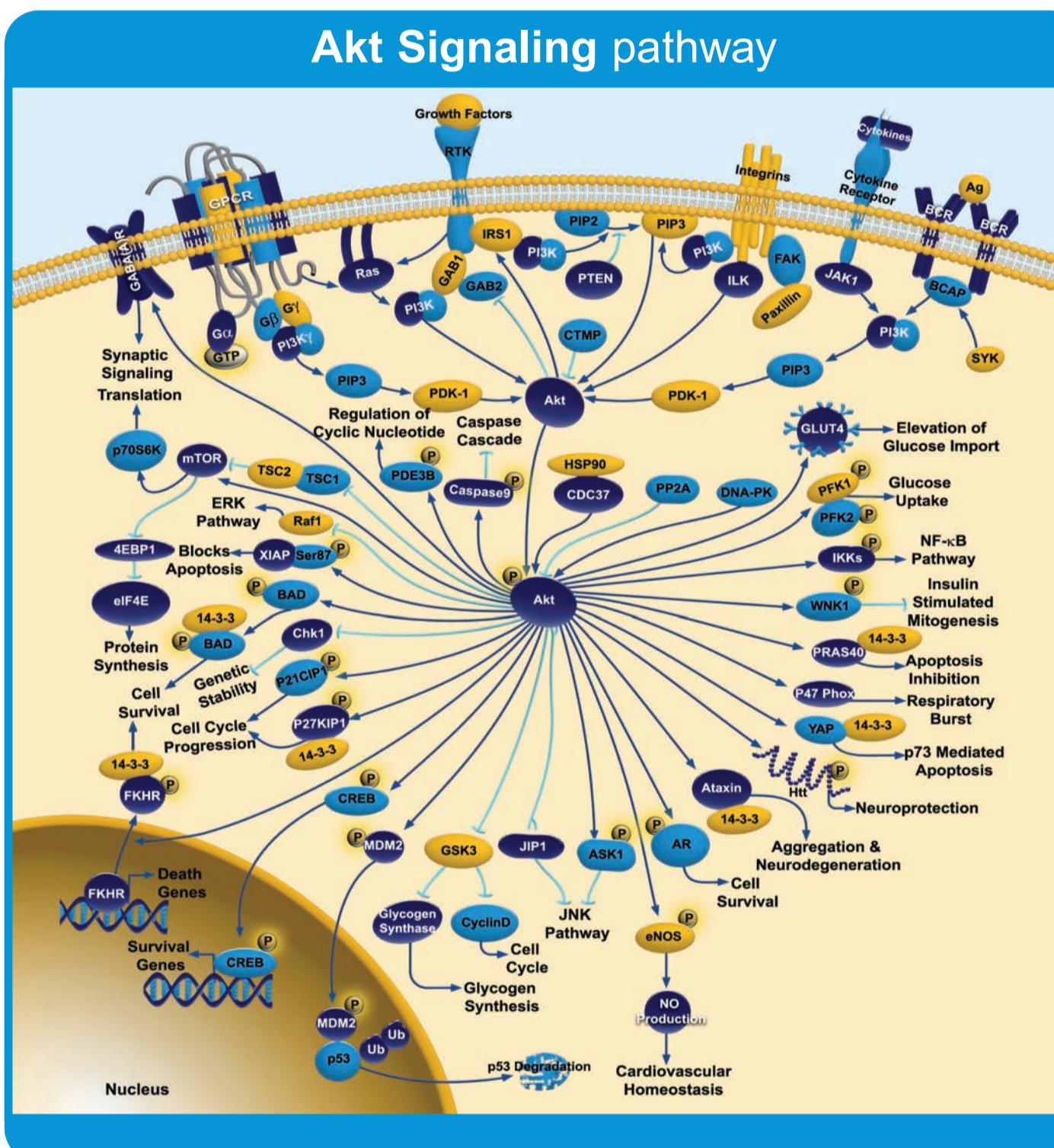
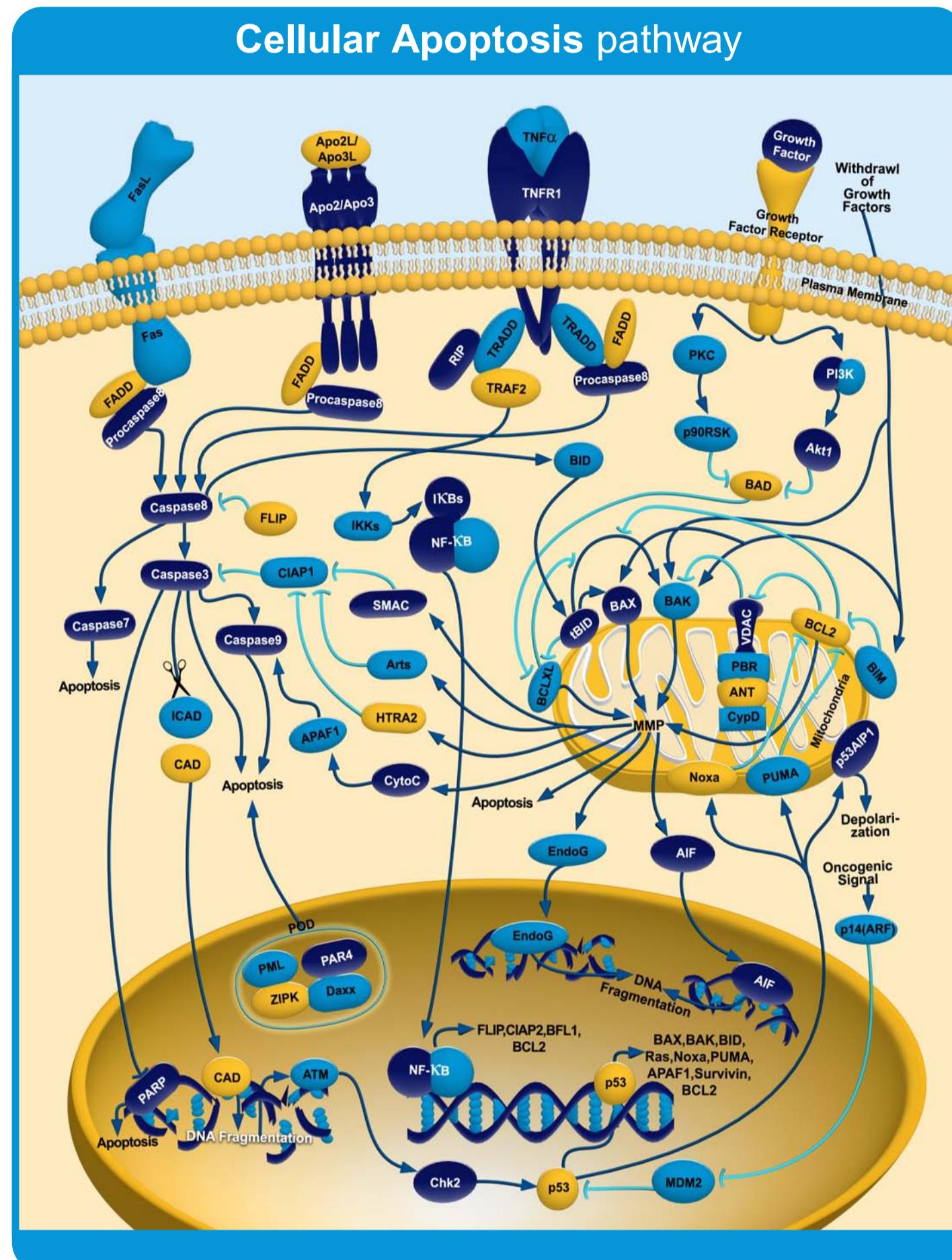


Cancer pathways



Product	Clonality	Applications tested	Host	Species reactivity	Datasheet
Active caspase 3	P	ICC, IHC-P IHC-Fo, Flow cyt	Rb	Hu	13847
ADAM12	P	WB	Rb	Hu	28225
ADAM17	P	WB	Rb	Hu, Mm, Rt	2051
AKT (phospho S473) [14-6]	M	ICC, WB	Rb	Hu, Mm	27773
AKT (phospho T308)	P	ELISA, IHC-P WB	Rb	Mm, Rt, Hu	38449
ATM (phospho S1981)	P	Flow cyt, ICC/IF, WB	Rb	Mm, Hu	2888
ATM [2C1 (1A1)]	M	IP, WB	Mm	Hu, Rt, Mm, Mk	78
Aurora A	P	ICC, IF, IHC-P WB	Rb	Hu	12875
Aurora B	P	ICC/IF, IF, IHC-P, WB	Rb	Hu, Mm, Rt	2254
B Raf [EP152Y]	M	Flow cyt, IHC-P, WB	Rb	Mm, Rt, Hu	33899
Bad [Y208]	P	IHC-P, WB	Rb	Hu, Rt, Mm	7977
Bax	P	IHC-P, IP, WB	Rb	Hu, Rt, Mm	7973
Bcl2	M	ICC, IP, WB	Mm	Hu, Mm, Rt, Rb	14389
Bmi1 [1.T.21]	P	IHC-P, HCPA, IP, WB	Rb	Hu, Rt, Mm	7963
c-Fos	P	ELISA, IHC-P, WB	Rb	Hu, Mm, Rt	31419
c-Jun	P	ChIP, IHC-Fr, IHC-P, WB	Mm	Hu, Mm, Ch	56
c-Myc [9E10] - ChIP Grade	M	WB	Rb	Hu, Mm, Rt	25758
Caspase 9	P	WB	Rb	Hu, Mm, Rt	21667
CCR2	P	WB	Rb	Hu	1416
E Cadherin [HECD-1]	M	IF, IHC-Fr, IHC-P, WB	Mm	Hu	11950
E2F1 [KH129]	M	WB	Mm	Hu	2430
EGFR	P	IHC-Fr, IHC-P, WB	Rb	Hu, Mm, D	2428
ErbB 2	P	IHC-Fr, IHC-P, WB	Rb	Hu, Rt, Mm	4819
ERK1 + ERK2 (phospho T185 + T202)	P	WB	Rb	Hu	4819
Estrogen Receptor alpha [33]	M	ICC, IHC-R, WB	Mm	Hu, Rt, Mm, C	2746
Estrogen Receptor beta	P	ICC/IF, IHC-Fr, IHC-P, WB	Rb	Hu, Mm, Rt, Sh	3577
Fibronectin [IST-9]	M	ICC, ICC/IF, IHC-Fr, IHC-P, WB	Mm	Hu, Mm, Ch, Pig, D, C	6328
HIF1 alpha [H1alpha67]	M	ICC/IF, IHC-Fr, IP, WB	Mm	Rt, Sh, Rb, C, Hu, Mm	1
HIF2 alpha	P	WB	Rb	Hu, Rt, Mm	199
Hsp27 [G3.1]	M	ELISA, IF, IHC-P, WB	Mm	Hu, Mm	2790
JNK1+NFK2 (phospho T183 + Y185)	P	IHC-Fr, WB	Rb	Hu, Mm	4821
MDM2 [2A10]	M	ICC, IHC-Fr, IHC-P, WB	Mm	Hu	16895
MMP2	P	IHC-Fr, IHC-P, WB	Rb	Mm, Rt, Ch, Hu	37150
MMP9	P	ELISA, IHC-Fr, IHC-P, WB	Rb	Mm, Hu	38898
mTOR [Y391]	M	Flow cyt, ICC/IF, IHC-P, WB	Rb	Hu, Mm, Rt	32028
MUC1 [C595 (NCRC48)]	M	Flow cyt, IHC-Fr, IHC-P, WB	Mm	Hu	28081
N Cadherin	P	ICC, IHC-Fr, IHC-P, WB	Rb	Mm, Rt, Ch, C, Hu	12221
NFKB p105 / p50 - ChIP Grade	P	ChIP, IHC-Fr, IHC-P, WB	Rb	Hu, Rt, Mm	7971
NFKB p65	P	ICC/IF, WB	Rb	Hu	16502
p19ARF	P	ELISA, ICC, ICC/IF, IHC-P, WB	Rb	Mm	80
p21	P	Flow cyt, ICC, IHC-Fr, WB	Rb	Hu, Rt, Mm	7960
p21 Ras [F132-62]	M	ICC, ICC/IF, IHC-Fr, IHC-P, WB	Mm	Hu, Mm, Rt	16907
p53 [Pab 240]	M	ICC, IHC-Fr, IHC-P, WB	Mm	Hu, Mm, Rt	26
pan Cadherin [mAbcam22744]	M	ICC/IF, WB	Mm	Hu, Rt	22744
PTEN [Y184]	M	Flow cyt, ICC, IHC-P, WB	Rb	Mm, Rt, Hu	32199
ROCK1 [EP786Y]	M	ICC/IF, IHC-P, WB	Rb	Mm, Rt, C, D, Hu	45171
Survivin	P	ICC/IF, WB	Rb	Hu, Mm	469
Telomerase	P	ICC/IF, WB	Rb	Hu, Mm	23699
TIE2 [C1_16]	M	ELISA, Flow cyt, WB	Mm	Hu	24859
TIMP1 [102D1]	M	IHC-P, IHC-Fo, WB	Mm	Hu	1827
TIMP2 - Loop 1	P	WB	Rb	Mm, Rt, C	38973
TNF alpha	P	ELISA, IHC-Fr, IHC-P, Neut, WB	Rb	Hu, Ch	6671
VEGF [VG-1]	M	ELISA, IHC-Fr, IHC-P, WB	Mm	Hu, Mm, Rt, D	1316
VEGF Receptor 1	P	IHC-Fr, IHC-P, WB	Rb	Hu, Rt, Mm	2350
VEGF Receptor 2	P	ICC/IF, IHC-Fr, IHC-P, WB	Rb	Hu, Mm, Rt, C, D	2349

www.abcam.com/cancer

Abcam in the USA

Abcam Inc.
1 Kendall Square, Ste 341
Cambridge,
MA 02139-1517 USA

Tel: 1-617-225-2272
Toll free: 1-888-77-ABCAM
Toll free Fax: 1-866-739-9884

Abcam in Europe

Abcam plc,
330 Cambridge Science Park
Milton Road,
Cambridge,
CB4 0FL UK

Tel: +44 1223 696000
Fax: +44 1223 771600

Abcam in Japan

Abcam K.K.
1-16-8 Nihonbashi Kakigaracho
Chuo-ku,
Tokyo 103-0014
Japan

Tel: +81-(0)3-6231-0940
Fax: +81-(0)3-6231-0941