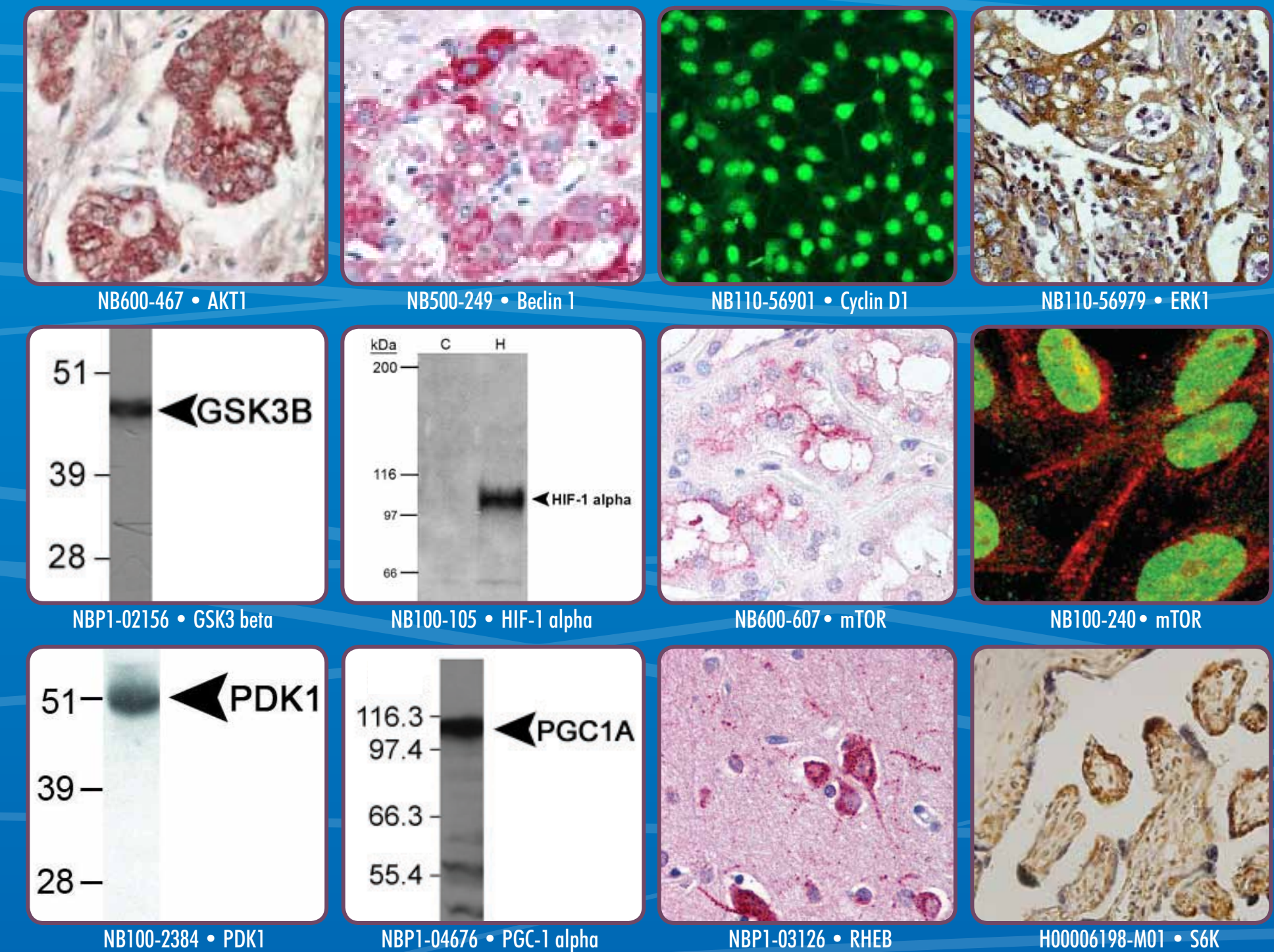


NOVUS
BIOLOGICALS®
Today's Research. Tomorrow's Discovery.



Catalog #	Product	Host	Clonality	Applications	Species
NB600-467	AKT1 Antibody	Rabbit	Polyclonal	ELISA, IF, IHC-P, WB	Hu
NB500-249	Bedin 1 Antibody	Rabbit	Polyclonal	ICC, IF, IHC-P, IP, WB	Hu, Mu
NB110-56901	Cyclin D1 (EP272Y) Antibody	Rabbit	Monoclonal	ICC, IHC, IP, WB	Hu, Rt
NB200-157	eIF4EBP1 Antibody	Rabbit	Polyclonal	WB	Hu, Mu, Po, Rt
NB110-56979	ERK1 Antibody (Y72)	Rabbit	Monoclonal	FACS, ICC, IHC-P, IP, WB	Hu, Mu, Rt
NBP1-02156	GSK3 beta Antibody	Rabbit	Polyclonal	WB	Hu, Mu, Rt, Sh, Ze
NB100-105	HIF-1 Alpha (H1alpha67)	Mouse	Monoclonal	ChIP, ICC, IF, IHC-Fr, IHC-P, IP, WB	Bv, Ft, Hu, Mk, Mu, Po, Rt, Sh, Rb
NB100-905	HIF-1 Alpha Antibody Pack	varies	varies	varies	Hu, Mu
NB600-607	mTOR [phospho Ser2448] Antibody	Rabbit	Polyclonal	ELISA, IHC-P, WB	Hu
NB100-240	mTOR Antibody	Rabbit	Polyclonal	ICC, IP	Hu, Rt
NB100-241	mTOR Antibody	Rabbit	Polyclonal	WB	Hu, Mu, Rt
NB100-932	mTOR Antibody Pack	Rabbit		IP, WB	Hu, Mu, Rt
NB100-2384	PDK1 Antibody	Rabbit	Polyclonal	WB	Bv, Ch, Hu, Mk, Mu, Po, Rt, Ze
NBP1-02583	PDK1 [phospho Ser241] Antibody	Rabbit	Polyclonal	ELISA, IHC-P	Hu
NBP1-04676	PGC-1 alpha Antibody	Rabbit	Polyclonal	WB	Ca, Eq, Hu, Mk, Mu, Rt
NB110-57347	PI3K Antibody (Y384)	Rabbit	Monoclonal	FACS, IP, WB	Hu
NB110-60524	PPAR gamma Antibody (8D1H8F4)	Mouse	Monoclonal	ELISA, IF, IHC-P, WB	Hu
NB100-2229	PTEN Antibody	Rabbit	Polyclonal	IP, WB	Hu, Mu
NB100-766	Raptor Antibody	Rabbit	Polyclonal	FACS, IP	Hu, Mu
NBP1-03126	RHEB Antibody	Rabbit	Polyclonal	IHC-P, WB	Hu, Mu, Rt
NB100-612	RICTOR Antibody	Rabbit	Polyclonal	FACS, IP, WB	Hu, Mu
H00006198-M01	S6K Antibody (1E12)	Mouse	Monoclonal	ELISA, IF, IHC-P, WB	Hu
NB110-40424	SIN1 Antibody	Rabbit	Polyclonal	IP, WB	Hu, Mu
NB100-2315	TSC1 Antibody	Rabbit	Polyclonal	IP, WB	Hu, Mu
NB110-57632	TSC2 Antibody (Y320)	Rabbit	Monoclonal	FACS, IHC, WB	Hu, Mu, Rt