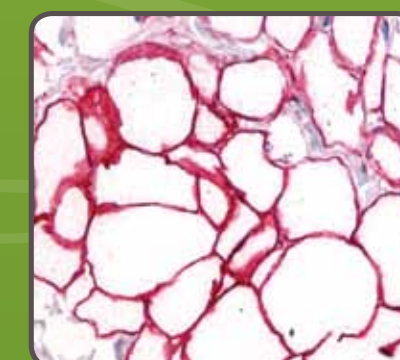
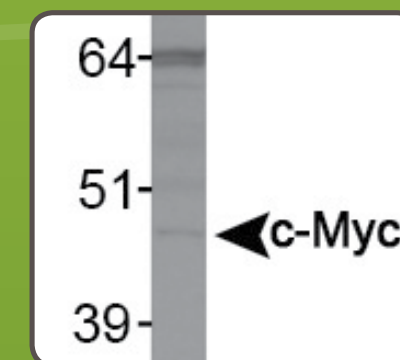
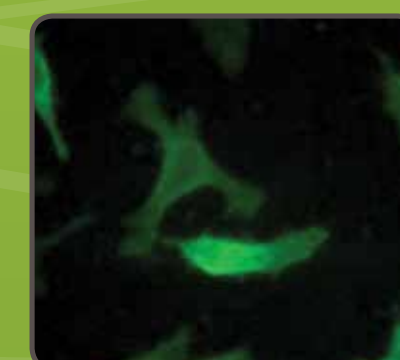
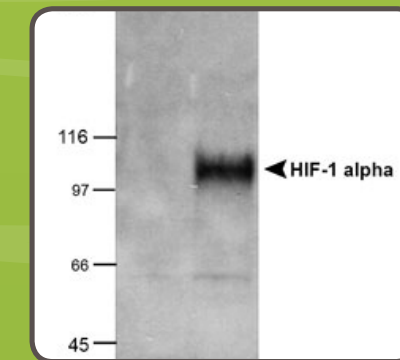
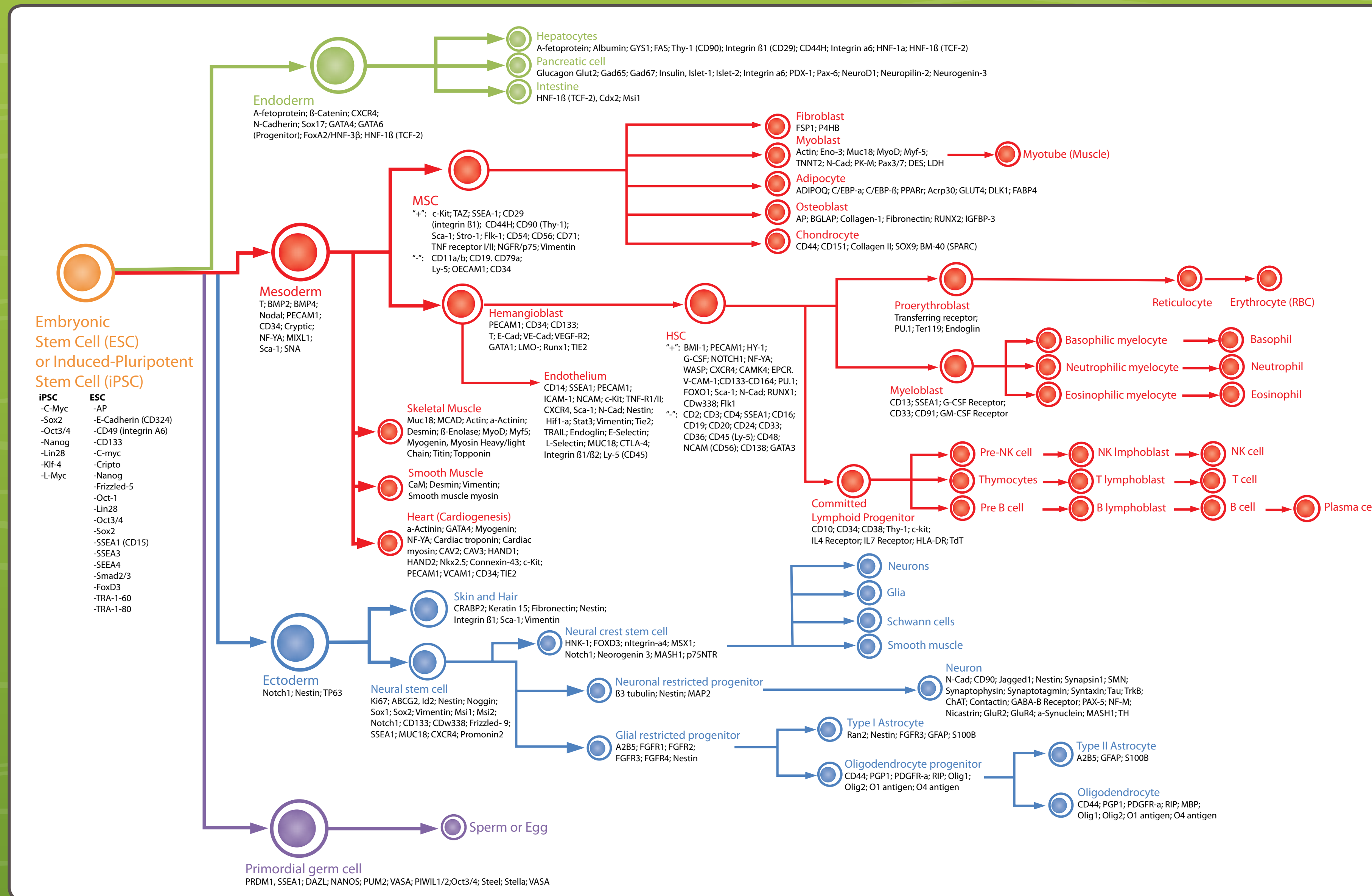


Stem Cell Differentiation Pathway



Antibodies for Stem Cell Differentiation Research

Product	Catalog#	Host	Clonality	Species	Applications
HIF-1 alpha Antibody (H1alpha67)	NB100-105	Mouse	Monoclonal	Bv, Fr, Hu, Mk, Mu, Po, Rb, Rt, Sh	ChIP, ICC, IF, IHC-Fr, IHC-P, IP, WB
HIF-1 beta Antibody (H1beta234)	NB100-124	Mouse	Monoclonal	Bv, Fr, Hu, Mu, Rt, Sh	IHC-P, WB
Tyrosine Hydroxylase Antibody	NB300-110	Sheep	Polyclonal	Hu, Mu, Rt	IF, IHC-Fr, WB
c-Myc Antibody (9E10)	NB600-302	Mouse	Monoclonal	Dr, Hu, Mu	ELISA, FACS, ICC, IF, IHC-Fr, IHC-P, IP, WB

Product	Catalog#	Host	Clonality	Species	Applications
c-Myc Antibody (9E11)	NB200-108	Mouse	Monoclonal	Ba, Ch, Hu, Mu	ChIP, ELISA, IHC-Fr, IHC-P, IP, WB
CD36 Antibody	NB400-144	Rabbit	Polyclonal	Bv, Hu, Mk, Mu, Rt	ICC, IF, IHC, WB
Synapsin I Antibody	NB300-104	Rabbit	Polyclonal	Hu, Mu, Rt	IF, IHC, IP, WB
SOX2 Antibody	NB110-37235	Rabbit	Polyclonal	Ch, Hu, Mu, Sh	FACS, ICC, IF, IHC-P, WB