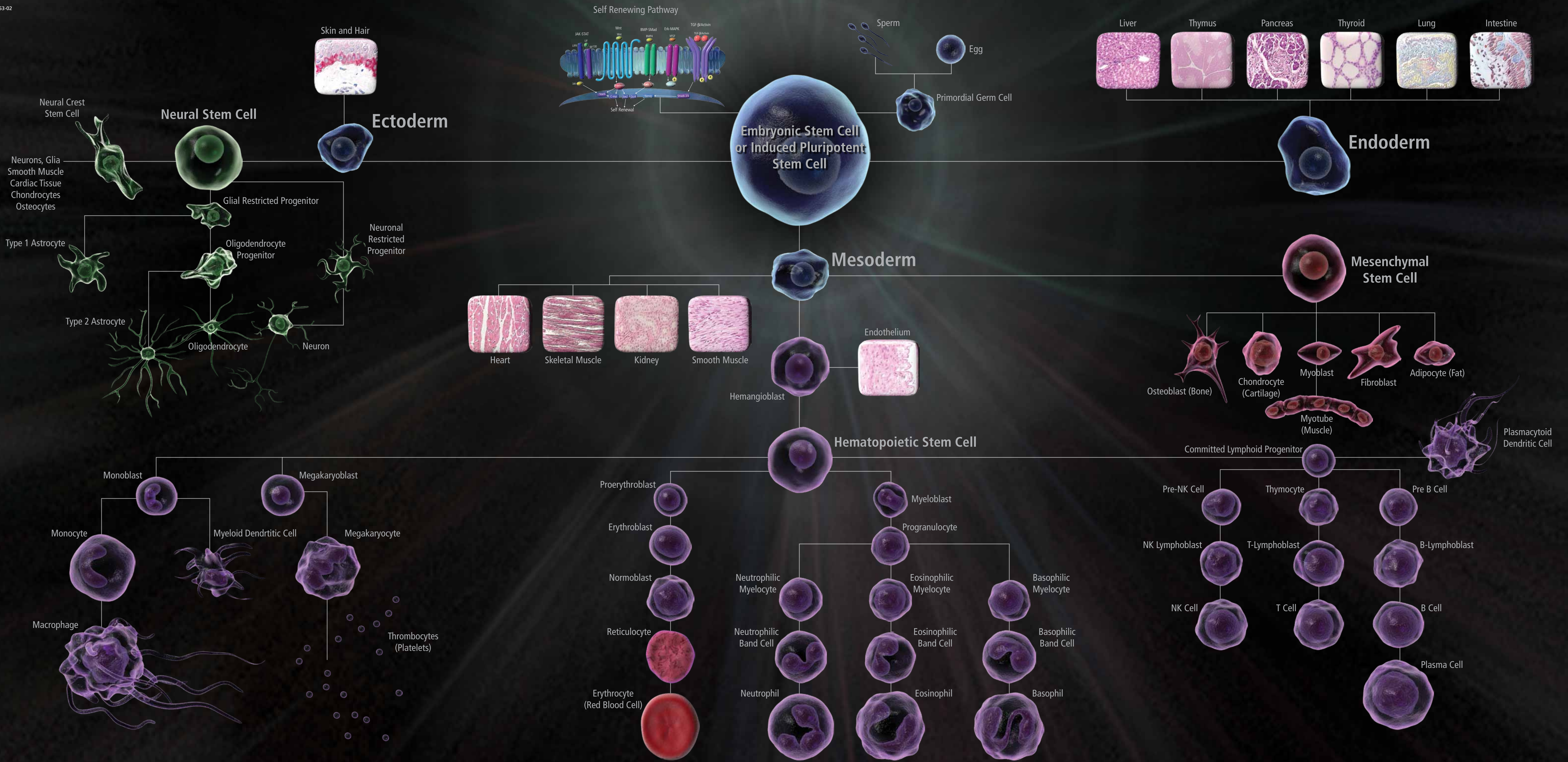


BD Stem Cell Resource

Markers of Self-Renewal and Differentiation



23-9953-02



White colored markers are available from BD Biosciences

Cell Type	Markers														
Embryonic Stem Cells	<table border="0"> <tr> <td>Neural Crest Cell Markers</td> <td> <ul style="list-style-type: none"> CD96 (Integrin α4) CD97 (Thy-1) CD98.1 (Thy-1.1) CD98.2 (Thy-1.2) </td> </tr> <tr> <td>Neural Stem Cell Markers</td> <td> <ul style="list-style-type: none"> CD133 (Prominin-1) CD134 (CXCR4) CD136 (VCAM-1) CD138 (ABCG2) CD146 (MCAM, MUC18) CD151 (Sialyl Lewis X) CD166 (NCAM, MUC18) CD184 (CXCR4) CD221 (p75, NGR/NTR) CD338 (ABCG2) </td> </tr> <tr> <td>Neuronal Markers</td> <td> <ul style="list-style-type: none"> MAP2B MASH1 β-tubulin (Tuj1) nestin NeuroD1 Neurofilament NF-M Neuron-restricted progenitors Neurogranin-2 Nicotin β5G1 Glial Pax-5 Pygycoprotein PSD-95 Semaphorin Receptor 5-HF 2AR SYN Syntaxin Synaptophysin Syntaxin Tau TfA8 Tubay Tyrosine Hydroxylase Vimentin </td> </tr> <tr> <td>Neuronal Stem Cell Markers</td> <td> <ul style="list-style-type: none"> CD133 (Prominin-1) CD134 (CXCR4) CD136 (VCAM-1) CD138 (ABCG2) CD146 (MCAM, MUC18) CD151 (Sialyl Lewis X) CD166 (NCAM, MUC18) CD184 (CXCR4) CD221 (p75, NGR/NTR) CD338 (ABCG2) </td> </tr> <tr> <td>Glial Restricted Progenitor</td> <td> <ul style="list-style-type: none"> CD133 (Prominin-1) CD134 (CXCR4) CD136 (VCAM-1) CD138 (ABCG2) CD146 (MCAM, MUC18) CD151 (Sialyl Lewis X) CD166 (NCAM, MUC18) CD184 (CXCR4) CD221 (p75, NGR/NTR) CD338 (ABCG2) </td> </tr> <tr> <td>Oligodendrocyte Markers</td> <td> <ul style="list-style-type: none"> CD117 (SCF R, c-kit) CD119 (TSP-1) CD133 (Prominin-1) CD134 (CXCR4) CD136 (VCAM-1) CD138 (ABCG2) CD146 (MCAM, MUC18) CD151 (Sialyl Lewis X) CD166 (NCAM, MUC18) CD184 (CXCR4) CD221 (p75, NGR/NTR) CD338 (ABCG2) </td> </tr> <tr> <td>Neuronal Restricted Progenitor</td> <td> <ul style="list-style-type: none"> CD133 (Prominin-1) CD134 (CXCR4) CD136 (VCAM-1) CD138 (ABCG2) CD146 (MCAM, MUC18) CD151 (Sialyl Lewis X) CD166 (NCAM, MUC18) CD184 (CXCR4) CD221 (p75, NGR/NTR) CD338 (ABCG2) </td> </tr> </table>	Neural Crest Cell Markers	<ul style="list-style-type: none"> CD96 (Integrin α4) CD97 (Thy-1) CD98.1 (Thy-1.1) CD98.2 (Thy-1.2) 	Neural Stem Cell Markers	<ul style="list-style-type: none"> CD133 (Prominin-1) CD134 (CXCR4) CD136 (VCAM-1) CD138 (ABCG2) CD146 (MCAM, MUC18) CD151 (Sialyl Lewis X) CD166 (NCAM, MUC18) CD184 (CXCR4) CD221 (p75, NGR/NTR) CD338 (ABCG2) 	Neuronal Markers	<ul style="list-style-type: none"> MAP2B MASH1 β-tubulin (Tuj1) nestin NeuroD1 Neurofilament NF-M Neuron-restricted progenitors Neurogranin-2 Nicotin β5G1 Glial Pax-5 Pygycoprotein PSD-95 Semaphorin Receptor 5-HF 2AR SYN Syntaxin Synaptophysin Syntaxin Tau TfA8 Tubay Tyrosine Hydroxylase Vimentin 	Neuronal Stem Cell Markers	<ul style="list-style-type: none"> CD133 (Prominin-1) CD134 (CXCR4) CD136 (VCAM-1) CD138 (ABCG2) CD146 (MCAM, MUC18) CD151 (Sialyl Lewis X) CD166 (NCAM, MUC18) CD184 (CXCR4) CD221 (p75, NGR/NTR) CD338 (ABCG2) 	Glial Restricted Progenitor	<ul style="list-style-type: none"> CD133 (Prominin-1) CD134 (CXCR4) CD136 (VCAM-1) CD138 (ABCG2) CD146 (MCAM, MUC18) CD151 (Sialyl Lewis X) CD166 (NCAM, MUC18) CD184 (CXCR4) CD221 (p75, NGR/NTR) CD338 (ABCG2) 	Oligodendrocyte Markers	<ul style="list-style-type: none"> CD117 (SCF R, c-kit) CD119 (TSP-1) CD133 (Prominin-1) CD134 (CXCR4) CD136 (VCAM-1) CD138 (ABCG2) CD146 (MCAM, MUC18) CD151 (Sialyl Lewis X) CD166 (NCAM, MUC18) CD184 (CXCR4) CD221 (p75, NGR/NTR) CD338 (ABCG2) 	Neuronal Restricted Progenitor	<ul style="list-style-type: none"> CD133 (Prominin-1) CD134 (CXCR4) CD136 (VCAM-1) CD138 (ABCG2) CD146 (MCAM, MUC18) CD151 (Sialyl Lewis X) CD166 (NCAM, MUC18) CD184 (CXCR4) CD221 (p75, NGR/NTR) CD338 (ABCG2)
Neural Crest Cell Markers	<ul style="list-style-type: none"> CD96 (Integrin α4) CD97 (Thy-1) CD98.1 (Thy-1.1) CD98.2 (Thy-1.2) 														
Neural Stem Cell Markers	<ul style="list-style-type: none"> CD133 (Prominin-1) CD134 (CXCR4) CD136 (VCAM-1) CD138 (ABCG2) CD146 (MCAM, MUC18) CD151 (Sialyl Lewis X) CD166 (NCAM, MUC18) CD184 (CXCR4) CD221 (p75, NGR/NTR) CD338 (ABCG2) 														
Neuronal Markers	<ul style="list-style-type: none"> MAP2B MASH1 β-tubulin (Tuj1) nestin NeuroD1 Neurofilament NF-M Neuron-restricted progenitors Neurogranin-2 Nicotin β5G1 Glial Pax-5 Pygycoprotein PSD-95 Semaphorin Receptor 5-HF 2AR SYN Syntaxin Synaptophysin Syntaxin Tau TfA8 Tubay Tyrosine Hydroxylase Vimentin 														
Neuronal Stem Cell Markers	<ul style="list-style-type: none"> CD133 (Prominin-1) CD134 (CXCR4) CD136 (VCAM-1) CD138 (ABCG2) CD146 (MCAM, MUC18) CD151 (Sialyl Lewis X) CD166 (NCAM, MUC18) CD184 (CXCR4) CD221 (p75, NGR/NTR) CD338 (ABCG2) 														
Glial Restricted Progenitor	<ul style="list-style-type: none"> CD133 (Prominin-1) CD134 (CXCR4) CD136 (VCAM-1) CD138 (ABCG2) CD146 (MCAM, MUC18) CD151 (Sialyl Lewis X) CD166 (NCAM, MUC18) CD184 (CXCR4) CD221 (p75, NGR/NTR) CD338 (ABCG2) 														
Oligodendrocyte Markers	<ul style="list-style-type: none"> CD117 (SCF R, c-kit) CD119 (TSP-1) CD133 (Prominin-1) CD134 (CXCR4) CD136 (VCAM-1) CD138 (ABCG2) CD146 (MCAM, MUC18) CD151 (Sialyl Lewis X) CD166 (NCAM, MUC18) CD184 (CXCR4) CD221 (p75, NGR/NTR) CD338 (ABCG2) 														
Neuronal Restricted Progenitor	<ul style="list-style-type: none"> CD133 (Prominin-1) CD134 (CXCR4) CD136 (VCAM-1) CD138 (ABCG2) CD146 (MCAM, MUC18) CD151 (Sialyl Lewis X) CD166 (NCAM, MUC18) CD184 (CXCR4) CD221 (p75, NGR/NTR) CD338 (ABCG2) 														