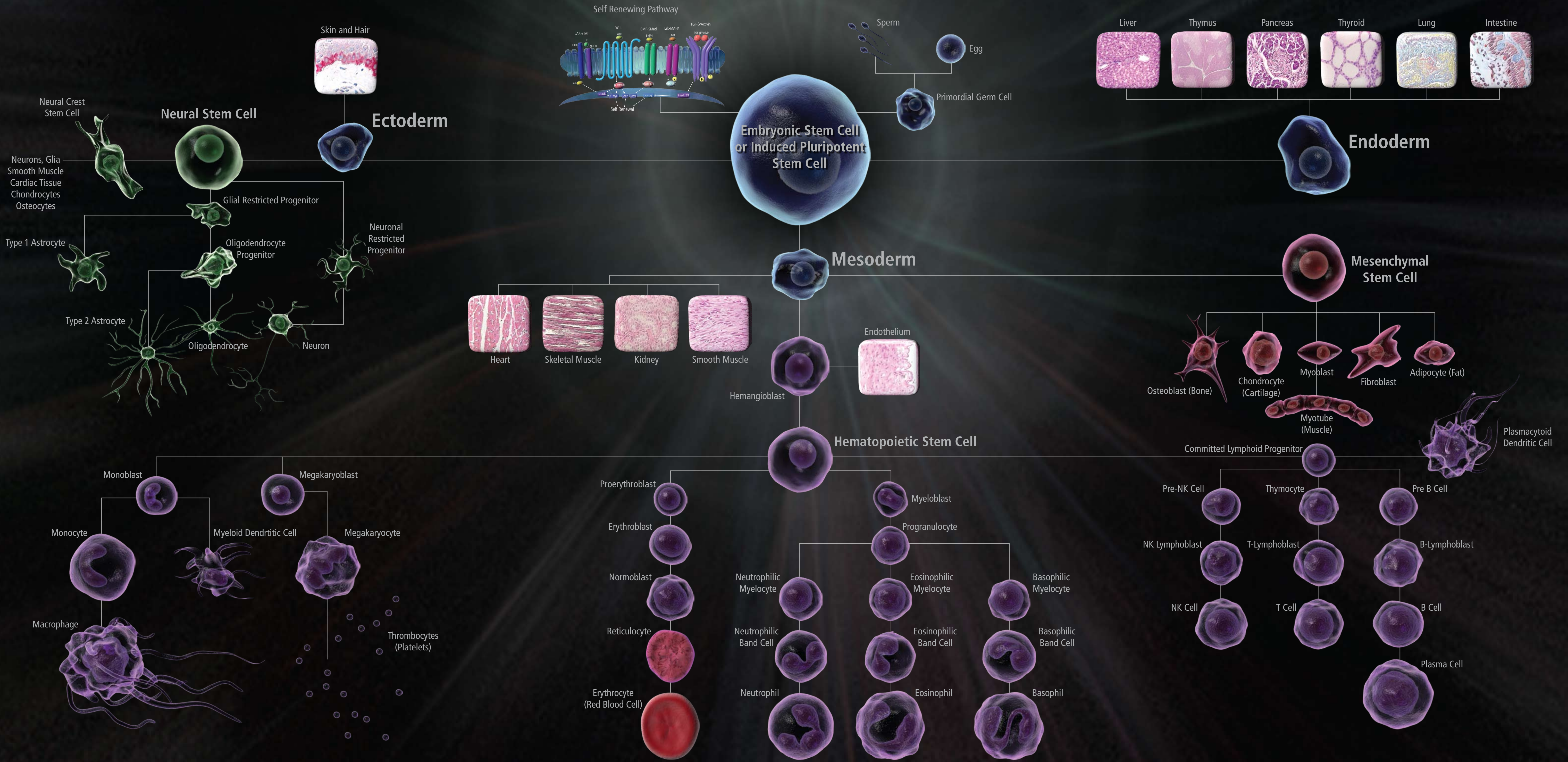


BD Stem Cell Resource

Markers of Self-Renewal and Differentiation



23-9953-01



| Ectoderm Markers | Embryonic Stem Cells | Mesoderm Markers | Endoderm Markers | Stem Cell Relevant Signaling |
|--|---|--|---|---|
| Astrocytes GFAP CD44 S100B GFAP3 Early Ectoderm Markers Nestin TGFβ Notch1 Neural Crest Cell Markers CD49b (Integrin α4) CD57 (HLK-1) CD271 (p75, NGR/NTR) Neural Stem Cell Markers CD15 (SSEA-1, Lewis X) CD155 (Stahl Lewis x) CD24 CD29 (Integrin β1) CD59 (Integrin α6) CD54 (ICAM-1) CD81 (TAPA-1) CD95 (Fas/PO-1) CD133 (AC133, Prominin-1) CD146 (MCAM, MUC18) CD184 (CXCR4) Sox1 CD271 (p75, NGR/NTR) CD349 (Riz2/Ed-9) | Neuronal Markers mGluR1 α-Synuclein (pY125) CD24 CD56 (NCAM) CD91 (TDP-1) CD90 (Thy-1) CD90.1 (Thy-1.1) CD92.2 (Thy-1.2) CD112 CD271 (p75, NGR/NTR) ChAT PSD-95 Doublecortin GABA B Receptor GAD65 GAP-43 (Neuromodulin) GluR2 GSK-3β GUS2/7 GUS2/7 Glutamine Synthetase Jagged1 MAP2 (α-hb) MAP2B Mash1 Oligodendrocyte Markers DT O4 CD146a (PDGFR α) MBP MBP Skin Precursor Cell Markers CRABP2 Fibronectin Nestin | Embryonic Stem Cell Markers Alkaline Phosphatase CD9 CD15 (SSEA-1, Lewis X) CD24 CD29 (Integrin β1) CD34 (Mucosalin, gp 105-120) CD30 CD49f (Integrin α6) CD133 (AC133, Prominin-1) CD194 (CXCR4) CD271 (p75, NGR/NTR) CD349 (Riz2/Ed-9) CD59 (Integrin α6) CD57 (HLK-1) CD271 (p75, NGR/NTR) Neuronal Markers mGluR1 α-Synuclein (pY125) CD24 CD56 (NCAM) CD91 (TDP-1) CD90 (Thy-1) CD90.1 (Thy-1.1) CD92.2 (Thy-1.2) CD112 CD271 (p75, NGR/NTR) ChAT PSD-95 Doublecortin GABA B Receptor GAD65 GAP-43 (Neuromodulin) GluR2 GSK-3β GUS2/7 GUS2/7 Glutamine Synthetase Jagged1 MAP2 (α-hb) MAP2B Mash1 Oligodendrocyte Markers DT O4 CD146a (PDGFR α) MBP MBP Skin Precursor Cell Markers CRABP2 Fibronectin Nestin | Early Mesoderm Markers BMP2 BMP4 CD31 (PECAM1) CD34 (Mucosalin, gp 105-120) CD49f CD49g CD49i CD49j CD49l CD49m CD49n CD49p CD49r CD49s CD49t CD49u CD49v CD49w CD49x CD49y CD49z CD49aa CD49ab CD49ac CD49ad CD49ae CD49af CD49ag CD49ah CD49ai CD49aj CD49ak CD49al CD49am CD49an CD49ao CD49ap CD49aq CD49ar CD49as CD49at CD49au CD49av CD49aw CD49ax CD49ay CD49az CD49ba CD49bb CD49bc CD49bd CD49be CD49bf CD49bg CD49bh CD49bi CD49bj CD49bk CD49bl CD49bm CD49bn CD49bo CD49bp CD49bq CD49br CD49bs CD49bt CD49bu CD49bv CD49bw CD49bx CD49by CD49bz CD49ca CD49cb CD49cc CD49cd CD49ce CD49cf CD49cg CD49ch CD49ci CD49cj CD49ck CD49cl CD49cm CD49cn CD49co CD49cp CD49cq CD49cr CD49cs CD49ct CD49cu CD49cv CD49cw CD49cx CD49cy CD49cz CD49da CD49db CD49dc CD49dd CD49de CD49df CD49dg CD49dh CD49di CD49dj CD49dk CD49dl CD49dm CD49dn CD49do CD49dp CD49dq CD49dr CD49ds CD49dt CD49du CD49dv CD49dw CD49dx CD49dy CD49dz CD49ea CD49eb CD49ec CD49ed CD49ee CD49ef CD49eg CD49eh CD49ei CD49ej CD49ek CD49el CD49em CD49en CD49eo CD49ep CD49eq CD49er CD49es CD49et CD49eu CD49ev CD49ew CD49ex CD49ey CD49ez CD49fa CD49fb CD49fc CD49fd CD49fe CD49ff CD49fg CD49fh CD49fi CD49fj CD49fk CD49fl CD49fm CD49fn CD49fo CD49fp CD49fq CD49fr CD49fs CD49ft CD49fu CD49fv CD49fw CD49fx CD49fy CD49fz CD49ga CD49gb CD49gc CD49gd CD49ge CD49gf CD49gg CD49gh CD49gi CD49gj CD49gk CD49gl CD49gm CD49gn CD49go CD49gp CD49gq CD49gr CD49gs CD49gt CD49gu CD49gv CD49gw CD49gx CD49gy CD49gz CD49ha CD49hb CD49hc CD49hd CD49he CD49hf CD49hg CD49hh CD49hi CD49hj CD49hk CD49hl CD49hm CD49hn CD49ho CD49hp CD49hq CD49hr CD49hs CD49ht CD49hu CD49hv CD49hw CD49hx CD49hy CD49hz CD49ia CD49ib CD49ic CD49id CD49ie CD49if CD49ig CD49ih CD49ii CD49ij CD49ik CD49il CD49im CD49in CD49io CD49ip CD49iq CD49ir CD49is CD49it CD49iu CD49iv CD49iw CD49ix CD49iy CD49iz CD49ja CD49jb CD49jc CD49jd CD49je CD49jf CD49jg CD49jh CD49ji CD49jj CD49jk CD49jl CD49jm CD49jn CD49jo CD49jp CD49jq CD49jr CD49js CD49jt CD49ju CD49jv CD49jw CD49jx CD49jy CD49jz CD49ka CD49kb CD49kc CD49kd CD49ke CD49kf CD49kg CD49kh CD49ki CD49kj CD49kl CD49km CD49kn CD49ko CD49kp CD49kq CD49kr CD49ks CD49kt CD49ku CD49kv CD49kw CD49kx CD49ky CD49kz CD49la CD49lb CD49lc CD49ld CD49le CD49lf CD49lg CD49lh CD49li CD49lj CD49lk CD49ll CD49lm CD49ln CD49lo CD49lp CD49lq CD49lr CD49ls CD49lt CD49lu CD49lv CD49lw CD49lx CD49ly CD49lz CD49ma CD49mb CD49mc CD49md CD49me CD49mf CD49mg CD49mh CD49mi CD49mj CD49mk CD49ml CD49mn CD49mo CD49mp CD49mq CD49mr CD49ms CD49mt CD49mu CD49mv CD49mw CD49mx CD49my CD49mz CD49na CD49nb CD49nc CD49nd CD49ne CD49nf CD49ng CD49nh CD49ni CD49nj CD49nk CD49nl CD49nm CD49nn CD49no CD49np CD49nq CD49nr CD49ns CD49nt CD49nu CD49nv CD49nw CD49nx CD49ny CD49nz CD49oa CD49ob CD49oc CD49od CD49oe CD49of CD49og CD49oh CD49oi CD49oj CD49ok CD49ol CD49om CD49on CD49oo CD49op CD49oq CD49or CD49os CD49ot CD49ou CD49ov CD49ow CD49ox CD49oy CD49oz CD49pa CD49pb CD49pc CD49pd CD49pe CD49pf CD49pg CD49ph CD49pi CD49pj CD49pk CD49pl CD49pm CD49pn CD49po CD49pp CD49pq CD49pr CD49ps CD49pt CD49pu CD49pv CD49pw CD49px CD49py CD49pz CD49qa CD49qb CD49qc CD49qd CD49qe CD49qf CD49qg CD49qh CD49qi CD49qj CD49qk CD49ql CD49qm CD49qn CD49qo CD49qp CD49qq CD49qr CD49qs CD49qt CD49qu CD49qv CD49qw CD49qx CD49qy CD49qz CD49ra CD49rb CD49rc CD49rd CD49re CD49rf CD49rg CD49rh CD49ri CD49rj CD49rk CD49rl CD49rm CD49rn CD49ro CD49rp CD49rq CD49rr CD49rs CD49rt CD49ru CD49rv CD49rw CD49rx CD49ry CD49rz CD49sa CD49sb CD49sc CD49sd CD49se CD49sf CD49sg CD49sh CD49si CD49sj CD49sk CD49sl CD49sm CD49sn CD49so CD49sp CD49sq CD49sr CD49ss CD49st CD49su CD49sv CD49sw CD49sx CD49sy CD49sz CD49ta CD49tb CD49tc CD49td CD49te CD49tf CD49tg CD49th CD49ti CD49tj CD49tk CD49tl CD49tm CD49tn CD49to CD49tp CD49tq CD49tr CD49ts CD49tt CD49tu CD49tv CD49tw CD49tx CD49ty CD49tz CD49ua CD49ub CD49uc CD49ud CD49ue CD49uf CD49ug CD49uh CD49ui CD49uj CD49uk CD49ul CD49um CD49un CD49uo CD49up CD49uq CD49ur CD49us CD49ut CD49uu CD49uv CD49uw CD49ux CD49uy CD49uz CD49va CD49vb CD49vc CD49vd CD49ve CD49vf CD49vg CD49vh CD49vi CD49vj CD49vk CD49vl CD49vm CD49vn CD49vo CD49vp CD49vq CD49vr CD49vs CD49vt CD49vu CD49vv CD49vw CD49vx CD49vy CD49vz CD49wa CD49wb CD49wc CD49wd CD49we CD49wf CD49wg CD49wh CD49wi CD49wj CD49wk CD49wl CD49wm CD49wn CD49wo CD49wp CD49wq CD49wr CD49ws CD49wt CD49wu CD49wv CD49ww CD49wx CD49wy CD49wz CD49xa CD49xb CD49xc CD49xd CD49xe CD49xf CD49xg CD49xh CD49xi CD49xj CD49xk CD49xl CD49xm CD49xn CD49xo CD49xp CD49xq CD49xr CD49xs CD49xt CD49xu CD49xv CD49xw CD49xx CD49xy CD49xz CD49ya CD49yb CD49yc CD49yd CD49ye CD49yf CD49yg CD49yh CD49yi CD49yj CD49yk CD49yl CD49ym CD49yn CD49yo CD49yp CD49yq CD49yr CD49ys CD49yt CD49yu CD49yv CD49yw CD49yx CD49yy CD49yz CD49za CD49zb CD49zc CD49zd CD49ze CD49zf CD49zg CD49zh CD49zi CD49zj CD49zk CD49zl CD49zm CD49zn CD49zo CD49zp CD49zq CD49zr CD49zs CD49zt CD49zu CD49zv CD49zw CD49zx CD49zy CD49zz | Definitive Endoderm Markers α-Fetoprotein B-Catenin CD184 (CXCR4) GATA4 GATA6 Hepatic Endoderm Markers α-Fetoprotein Albumin CD29 (Integrin β1) CD44f (Pgp-1, ICAM) CD49f (Integrin α6) Pancreatic Endoderm Markers CD49f (Integrin α6) CD49g CD49h CD49i CD49j CD49k CD49l CD49m CD49n CD49o CD49p CD49q CD49r CD49s CD49t CD49u CD49v CD49w CD49x CD49y CD49z CD49aa CD49ab CD49ac CD49ad CD49ae CD49af CD49ag CD49ah CD49ai CD49aj CD49ak CD49al CD49am CD49an CD49ao CD49ap CD49aq CD49ar CD49as CD49at CD49au CD49av CD49aw CD49ax CD49ay CD49az CD49ba CD49bb CD49bc CD49bd CD49be CD49bf CD49bg CD49bh CD49bi CD49bj CD49bk CD49bl CD49bm CD49bn CD49bo CD49bp CD49bq CD49br CD49bs CD49bt CD49bu CD49bv CD49bw CD49bx CD49by CD49bz CD49ca CD49cb CD49cc CD49cd CD49ce CD49cf CD49cg CD49ch CD49ci CD49cj CD49ck CD49cl CD49cm CD49cn CD49co CD49cp CD49cq CD49cr CD49cs CD49ct CD49cu CD49cv CD49cw CD49cx CD49cy CD49cz CD49da CD49db CD49dc CD49dd CD49de CD49df CD49dg CD49dh CD49di CD49dj CD49dk CD49dl CD49dm CD49dn CD49do CD49dp CD49dq CD49dr CD49ds CD49dt CD49du CD49dv CD49dw CD49dx CD49dy CD49dz CD49ea CD49eb CD49ec CD49ed CD49ee CD49ef CD49eg CD49eh CD49ei CD49ej CD49ek CD49el CD49em CD49en CD49eo CD49ep CD49eq CD49er CD49es CD49et CD49eu CD49ev CD49ew CD49ex CD49ey CD49ez CD49fa CD49fb CD49fc CD49fd CD49fe CD49ff CD49fg CD49fh CD49fi CD49fj CD49fk CD49fl CD49fm CD49fn CD49fo CD49fp CD49fq CD49fr CD49fs CD49ft CD49fu CD49fv CD49fw CD49fx CD49fy CD49fz CD49ga CD49gb CD49gc CD49gd CD49ge CD49gf CD49gg CD49gh CD49gi CD49gj CD49gk CD49gl CD49gm CD49gn CD49go CD49gp CD49gq CD49gr CD49gs CD49gt CD49gu CD49gv CD49gw CD49gx CD49gy CD49gz CD49ha CD49hb CD49hc CD49hd CD49he CD49hf CD49hg CD49hh CD49hi CD49hj CD49hk CD49hl CD49hm CD49hn CD49ho CD49hp CD49hq CD49hr CD49hs CD49ht CD49hu CD49hv CD49hw CD49hx CD49hy CD49hz CD49ia CD49ib CD49ic CD49id CD49ie CD49if CD49ig CD49ih CD49ii CD49ij CD49ik CD49il CD49im CD49in CD49io CD49ip CD49iq CD49ir CD49is CD49it CD49iu CD49iv CD49iw CD49ix CD49iy CD49iz CD49ja CD49jb CD49jc CD49jd CD49je CD49jf CD49jg CD49jh CD49ji CD49jj CD49jk CD49jl CD49jm CD49jn CD49jo CD49jp CD49jq CD49jr CD49js CD49jt CD49ju CD49jv CD49jw CD49jx CD49jy CD49jz CD49ka CD49kb CD49kc CD49kd CD49ke CD49kf CD49kg CD49kh CD49ki CD49kj CD49kl CD49km CD49kn CD49ko CD49kp CD49kq CD49kr CD49ks CD49kt CD49ku CD49kv CD49kw CD49kx CD49ky CD49kz CD49la CD49lb CD49lc CD49ld CD49le CD49lf CD49lg CD49lh CD49li CD49lj CD49lk CD49ll CD49lm CD49ln CD49lo CD49lp CD49lq CD49lr CD49ls CD49lt CD49lu CD49lv CD49lw CD49lx CD49ly CD49lz CD49ma CD49mb CD49mc CD49md CD49me CD49mf CD49mg CD49mh CD49mi CD49mj CD49mk CD49ml CD49mn CD49mo CD49mp CD49mq CD49mr CD49ms CD49mt CD49mu CD49mv CD49mw CD49mx CD49my CD49mz CD49na CD49nb CD49nc CD49nd CD49ne CD49nf CD49ng CD49nh CD49ni CD49nj CD49nk CD49nl CD49nm CD49nn CD49no CD49np CD49nq CD49nr CD49ns CD49nt CD49nu CD49nv CD49nw CD49nx CD49ny CD49nz CD49oa CD49ob CD49oc CD49od CD49oe CD49of CD49og CD49oh CD49oi CD49oj CD49ok CD49ol CD49om CD49on CD49oo CD49op CD49oq CD49or CD49os CD49ot CD49ou CD49ov CD49ow CD49ox CD49oy CD49oz CD49pa CD49pb CD49pc CD49pd CD49pe CD49pf CD49pg CD49ph CD49pi CD49pj CD49pk CD49pl CD49pm CD49pn CD49po CD49pp CD49pq CD49pr CD49ps CD49pt CD49pu CD49pv CD49pw CD49px CD49py CD49pz CD49qa CD49qb CD49qc CD49qd CD49qe CD49qf CD49qg CD49qh CD49qi CD49qj CD49qk CD49ql CD49qm CD49qn CD49qo CD49qp CD49qq CD49qr CD49qs CD49qt CD49qu CD49qv CD49qw CD49qx CD49qy CD49qz CD49ra CD49rb CD49rc CD49rd CD49re CD49rf CD49rg CD49rh CD49ri CD49rj CD49rk CD49rl CD49rm CD49rn CD49ro CD49rp CD49rq CD49rr CD49rs CD49rt CD49ru CD49rv CD49rw CD49rx CD49ry CD49rz CD49sa CD49sb CD49sc CD49sd CD49se CD49sf CD49sg CD49sh CD49si CD49sj CD49sk CD49sl CD49sm CD49sn CD49so CD49sp CD49sq CD49sr CD49ss CD49st CD49su CD49sv CD49sw CD49sx CD49sy CD49sz CD49ta CD49tb CD49tc CD49td CD49te CD49tf CD49tg CD49th CD49ti CD49tj CD49tk CD49tl CD49tm CD49tn CD49to CD49tp CD49tq CD49tr CD49ts CD49tt CD49tu CD49tv CD49tw CD49tx CD49ty CD49tz CD49ua CD49ub CD49uc CD49ud CD49ue CD49uf CD49ug CD49uh CD49ui CD49uj CD49uk CD49ul CD49um CD49un CD49uo CD49up CD49uq CD49ur CD49us CD49ut CD49uu CD49uv CD49uw CD49ux CD49uy CD49uz CD49va CD49vb CD49vc CD49vd CD49ve CD49vf CD49vg CD49vh CD49vi CD49vj CD49vk CD49vl CD49vm CD49vn CD49vo CD49vp CD49vq CD49vr CD49vs CD49vt CD49vu CD49vv CD49vw CD49vx |