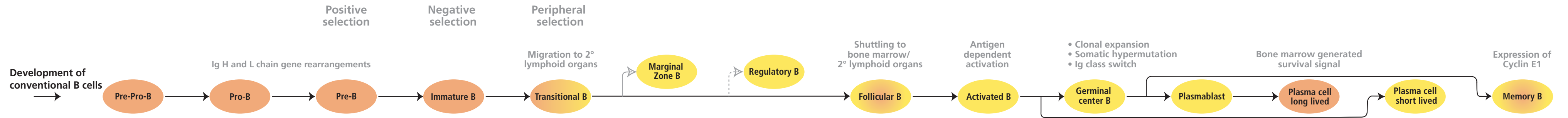
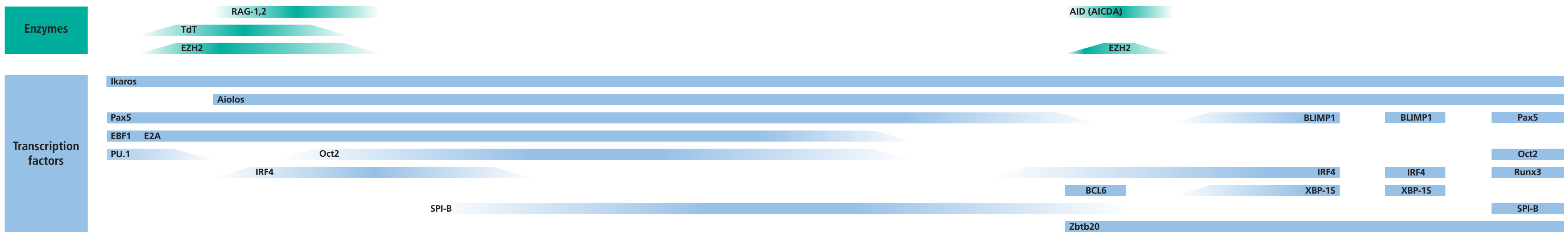


23-17817-00



Immunoglobulin	IgM ⁻	IgM ⁻	IgM ⁻	IgM ⁺ , IgD ⁻	IgM ^{high} , IgD ^{low}	IgM ⁺ , IgD ^{low}	IgM ^{high} , IgD ^{var}	IgM ^{low} , IgD ⁺	IgM ⁺ , IgD ⁺	IgM/G/A/E ⁺ , IgD ^{var}	IgM/G/A/E ⁺	Ig ⁻	Ig ⁻	IgM/G/A/E ⁺ , IgD ⁻
Mouse positive markers	CD43, CD93, B220, CXCR4, Flt3, IL7R	CD19, CD24, CD43, B220, IL7R	CD19, CD24, CD25 ^{var} , B220, BP1, IL7R, Siglec-G	CD19, CD24, CD93, CD268, B220	CD19, CD21 ^{var} , CD23 ^{var} , CD24, CD93, CD268, B220, CCR7	CD1d, CD9, CD21 ^{high} , CD22 ^{high} , CD35 ^{high} , B220	CD1d ^{high} , CD5, CD19, CD24, TIM-1 [±]	CD19, CD22, CD23, CD38, CD268, B220, CCR7, CXCR5	CD27, CD69, CD80, CD95, CD268, B220, CXCR5, Flt3, MHCII ^{high}	CD19, CD20, CD37, CD95, CD268, CXCR5, GL7, Siglec2	CD19, CD138, CD268, CXCR4, MHCII	CD28, CD138, CD268, CD269, CXCR4 ^{high}	CD28, CD93, CD138, CD268, CD269, CXCR4 ^{high}	CD38 ^{var} , CD62L ^{var} , CD80 ^{var} , CD95 ^{low} , CD268, B220, CCR7, CXCR5
Mouse negative markers	CD19, CD24 ^{low} , ckit ^{low}	BP1, ckit ^{low} , Flt3	CD43 ^{low} , ckit	CD23, CD43	CD23, CD93	CD23, CD93	CD62L, CD93 ^{var}	CD1d ^{low} , CD21/35 ^{low} , CD93	CD138, CXCR4	CD38 ^{low} , CD93	B220 ^{low} , Flt3	CD19, CD38 ^{low} , B220 ^{low} , MHCII	CD19, CD38 ^{low} , B220 ^{low} , MHCII ^{low}	
Human positive markers	CD10, CD34, CD38	CD10, CD19, CD24, CD34, CD38, IL7/3R	CD10, CD19, CD20, CD24, CD38, IL7/4/3R	CD10, CD19, CD20, CD21, CD24 ^{high} , CD38 ^{high} , CD40, CD268, IL4R	CD5, CD19, CD20, CD21, CD24 ^{high} , CD38 ^{high} , CD268, CCR7	CD1c, CD19, CD20, CD21 ^{high} , CD27 ^{var}	CD1d ^{high} , CD5, CD19, CD21, CD24 ^{high}	CD19, CD20, CD21, CD22, CD23, CD24, CD268, CCR7, CXCR5	CD19, CD20, CD25, CD27, CD30, CD69, CD80, CD86, CD95, CD135, CD268, CXCR5	CD10, CD19, CD20, CD23, CD27, CD38 ^{high} , CD95, CD268, CD269, CXCR5	CD19, CD27 ^{high} , CD38 ^{high} , CD268, CD269, MHCII	CD27, CD28, CD38 ^{high} , CD138, CD268, CD269, CXCR4	CD27, CD28, CD38 ^{high} , CD138, CD268, CD269, CXCR4	CD19, CD20, CD27 ^{var} , CD40, CD268, CCR7, CXCR4,5,7
Human negative markers	ckit ^{low}	ckit ^{low}	CD34, ckit	CD27, ckit, IL7R	CD10 ^{low} , CD27		CD27 ^{var}	CD10, CD27, CD24 ^{low} , CD38 ^{low}		CD24 ^{low}	CD20, CD138	CD19 ^{low} , CD20, MHCII	CD19 ^{low} , CD20, MHCII ^{low}	CD23 ^{low} , CD38



Bone marrow location (orange circle) 2° lymphoid organ location (yellow circle) Cells circulating in both marrow and lymphoid locations (yellow-orange circle)