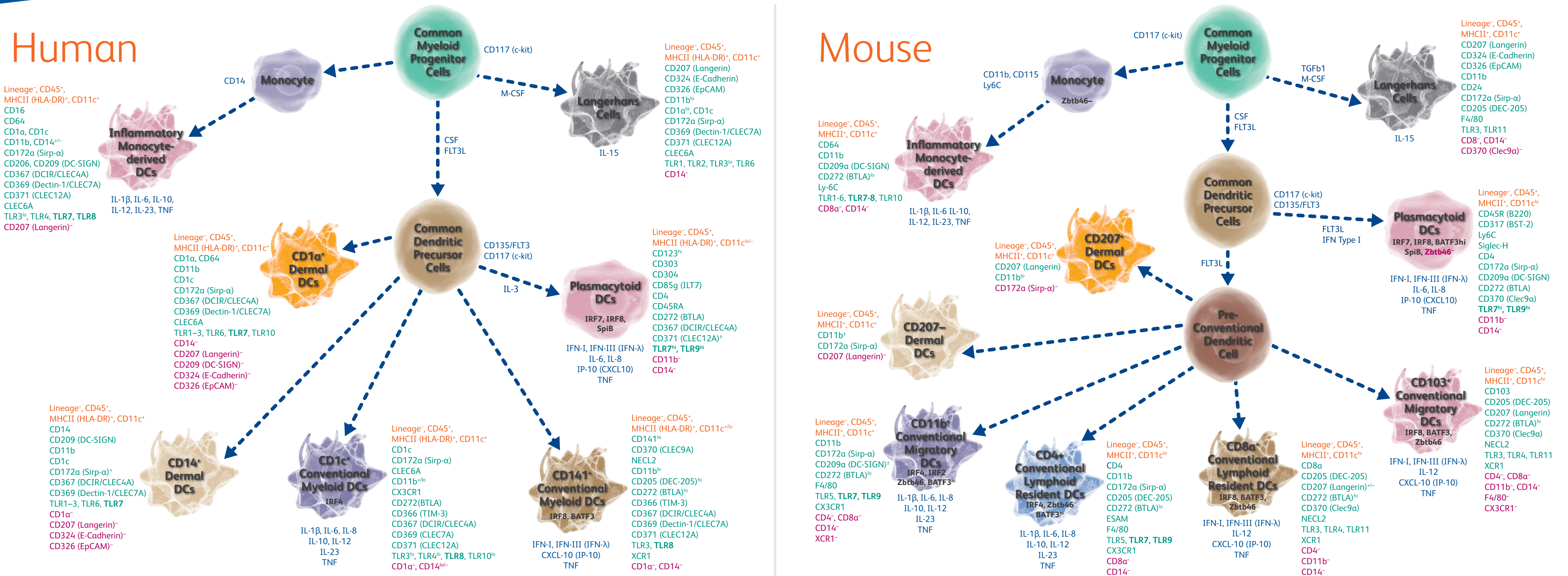


Overview of Dendritic Cell Development



For Research Use Only. Not for use in diagnostic or therapeutic procedures.
23-19088-00

BD Life Sciences, San Jose, CA, 95131, USA

bdbiosciences.com

© 2016 BD. BD and the BD Logo are trademarks of Becton, Dickinson and Company.

KEY

Core dendritic cell phenotyping markers
Lineage (Hu): CD3⁻, CD19⁻, CD20⁻, CD56⁻, CD14⁺
Lineage (Ms): CD3⁻, CD19⁻, CD49b⁻ or NK1.1⁻, CD14⁺

Dendritic cell subset markers (positive)
TLR7, TLR8 and TLR9 are endosomal and require intracellular staining

Dendritic cell subset markers (negative)

Secreted cytokines

Transcription factors

Markers on this schematic can be altered as a result of cellular environment, differentiation state and other factors. DC biomarker profiles were obtained from RNA and protein level analysis provided in peer reviewed publications and public databases. Visit bdbiosciences.com/DC for more information.

¹ CD14 is negative or low in all Human DC subsets except for CD14⁺ Dermal DCs and inflammatory DCs.

