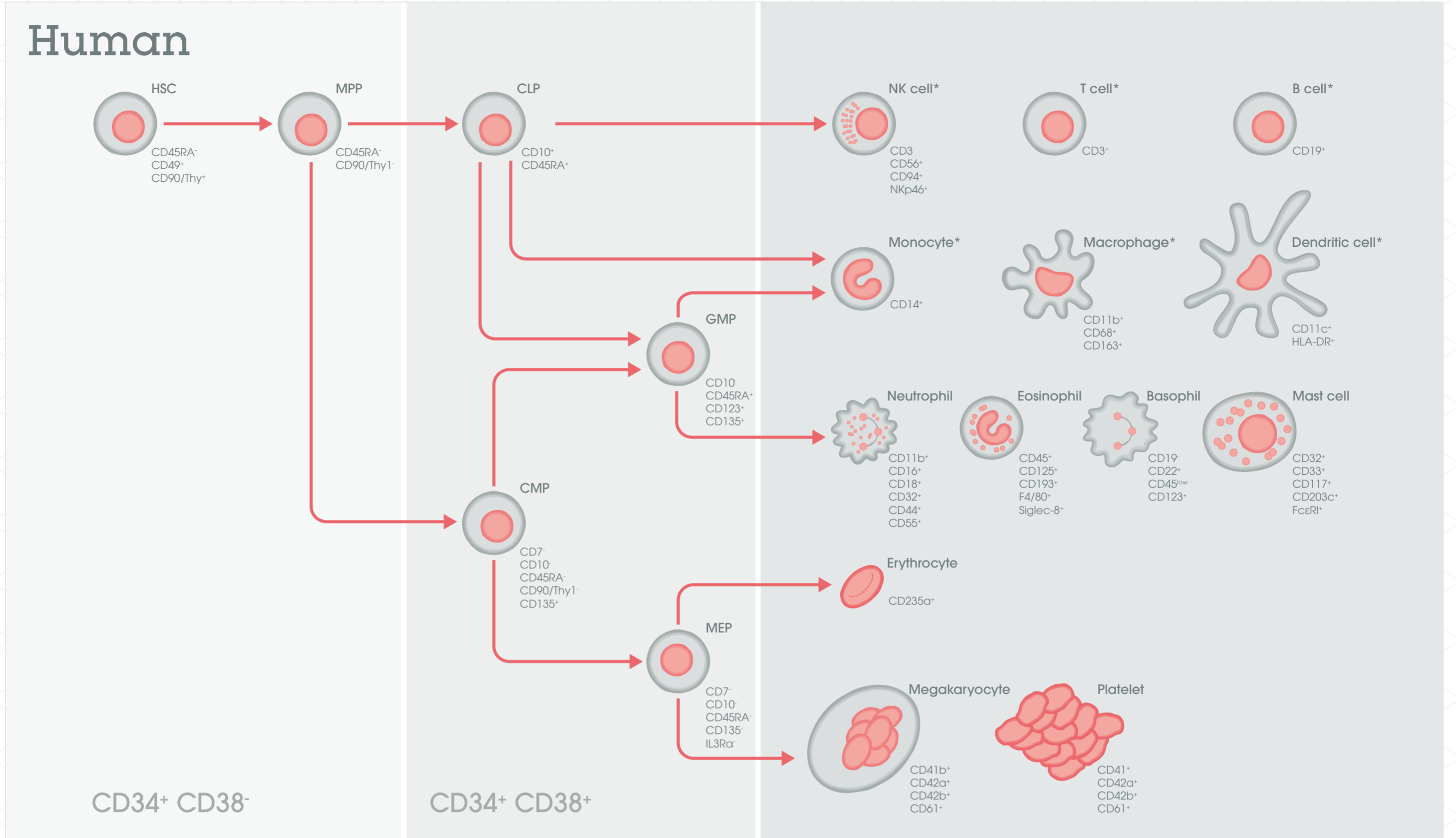


Immune cell markers

abcam

+



+

