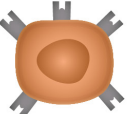


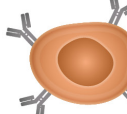
Immune Cell Phenotyping Key Markers

Lymphocytes



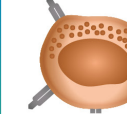
T Cell

- CD3
- CD4
- CD8



B Cell

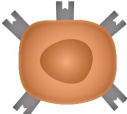
- CD19
- CD20
- CD21



Natural Killer Cell

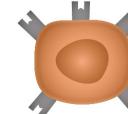
- CD56
- CD335
- CD8

CD3⁺ CD4⁺ T Cells



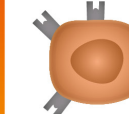
Naïve

- CD45RA
- CCR7
- CD62L



Regulatory

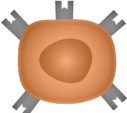
- FOXP3
- CD103



Memory

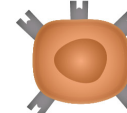
- CD25
- CD45RO
- CD62L

CD3⁺ CD8⁺ T Cells



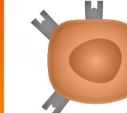
Naïve

- CD45RA



Cytotoxic

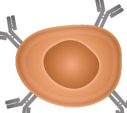
- Granzyme A
- Granzyme B



Memory

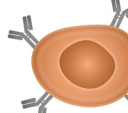
- CD45RO

CD19⁺ CD20⁺ B cells



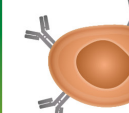
Follicular

- CD21
- CD22
- CD23



Activated

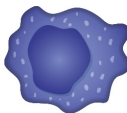
- CD25
- CD27
- CD69



Memory

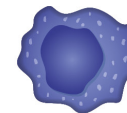
- CD40
- CXCR4
- CXCR5

CD14⁺ CD11b⁺ Macrophages



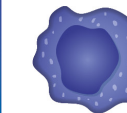
M1

- CD68
- CD86
- CD282



M2

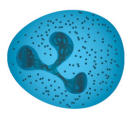
- CD163
- CD200R
- CD206



TAM

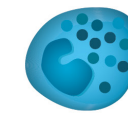
- CD81
- CD106
- Dectin-1

Granulocytes



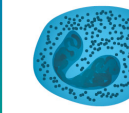
Neutrophils

- CD13
- CD15
- CD33



Basophils

- CD69
- CD123
- Fcε R1α



Eosinophils

- CD9
- CD23
- Siglec 8

Apoptosis Analysis by Western Blotting

Intrinsic Pathway

Early

- Bcl-2
- BID
- BAX
- PUMA

Mid

- Cytochrome C
- APAF-1
- Caspase-9

Late

- DFF40
- DFF45
- PARP-1

Extrinsic Pathway

Early

- RIPK3
- RIPK4
- FADD
- Caspase-8

Mid

- Caspase-3
- Caspase-7

Late

- DFF40
- DFF45
- PARP-1

Apoptosis Analysis by Flow Cytometry

Intrinsic Pathway

Early

- Bcl-2

Mid

- Mitochondrial Membrane Potential Kit
- Mitochondrial Permeability Transition Kit

Late

- DAPI
- Hoechst
- PI
- 7AAD

Extrinsic Pathway

Early

- Caspase-3 /7 Kits
- Caspase-8 Kits

Mid

- Caspase-9*
- Poly Caspase*
- Annexin V*
- FLISP Serine Protease*
- Cathepsin B*

*Kits

Late

- DAPI
- Hoechst
- PI
- 7AAD