免疫学



ヒト末梢単核細胞中の主要な白血球細胞 サブセットの詳細な特性解析

45 色フルスペクトル Agilent NovoCyte Opteon スペクトル フローサイトメータの利用

著者

Ming Lei, Laurissa Ouaguia, Garret Guenther, Peifang Ye, Nancy Li, Agilent Technologies, Inc.

概要

フローサイトメトリーは、免疫細胞のさまざまな細胞サブポピュレーションを詳細に分析し、免疫システムの包括的な評価を行う強力な手法です。スペクトルフローサイトメトリーにより性能が向上し、より多くのマーカー、より高い柔軟性、より優れたデータ分解能をサポートできるようになりました。このアプリケーションノートでは、末梢血単核細胞(PBMC)サンプル内の免疫細胞の分布を分析する Agilent NovoCyte Opteon スペクトルフローサイトメータ用に設計された 45 色スペクトルフローサイトメトリーパネルについてご紹介します。

はじめに

免疫システムは、非常に複雑な細胞ネットワークを用いて、死んだ細胞 や不完全な細胞を認識して除去すると同時に、病原体から身体を守るこ とができます。末梢血細胞は、免疫システムの主要な促進因子として、活 性化、増殖、および一連の特定のマーカー発現と働きによって特徴付け られるさまざまなサブセットへの分化を起こします。末梢血は収集が簡単 で非侵襲性であるため、末梢血単核細胞(PBMC)の供給源として広く 入手可能で、基礎研究、トランスレーショナルリサーチ、臨床応用で広く 使用されています¹³。健康なドナーの場合、PBMC は、リンパ球(T細胞、 B細胞、ナチュラルキラー (NK)細胞、NKT細胞、γδT細胞、自然リン パ球(ILC)を含む)、単球、好塩基球、樹状細胞(DC)など、いくつか の免疫細胞タイプで構成された異種細胞集団となります^{1,3,4}。

末梢血中のリンパ球の多くは、CD3+を発現する T 細胞(PBMC の 45 ~ 70 %) で構成されています。 $\alpha\beta$ TCR または $\gamma\delta$ TCR の発現に基づい て、2 種類の T 細胞を区別できます⁵。αβTCR を発現する T 細胞は、さ らに従来の CD4+ T ヘルパー細胞(25~60%)と CD8+ T 細胞傷害 性細胞(5~30%)に細分化できます。健康なドナー PBMC の CD4+ T 細胞とCD8+ T 細胞の比率は約 2:1 です^{3,6}。CD4+ T ヘルパーは免 疫恒常性と炎症の重要なメディエーターです。特定の表面マーカー /タ ンパク質、サイトカイン、ケモカイン受容体、特定の転写因子の発現プ ロファイルに基づいて、CD4+T細胞はさらにTh1、Th2、Th17、Th9、 Th22、濾胞ヘルパー(Tfh)細胞、胸腺から新たに移出された T 細胞 (RTE)、および制御性 T 細胞 (Treg) サブセットに分類されます⁷⁻¹⁰。 さらに、抗原提示細胞は、ナイーブ T細胞に抗原を提示した後に活性化 され、エフェクター T 細胞とメモリー細胞に分化します。エフェクター細 胞は感染部位に移動して病原体を排除します。これらのエフェクター細 胞は短命ですが、形成されたメモリー細胞サブセットは長期間生存する 可能性があります^{7,11}。従来の CD8+ T 細胞傷害性細胞は、細胞内感染 や腫瘍細胞と戦う適応免疫システムの重要な構成要素です。CD4+ T 細 胞と同様に、ナイーブ CD8+ T 細胞の活性化は、増殖と分化に関するほ ぼ自律的なプログラムを誘導し、エフェクター細胞とメモリー細胞の両 方を生成します⁷。免疫記憶は感染からの長期的な免疫と保護に非常に 重要であるため、メモリー T 細胞は免疫の重要な促進因子です⁷。相互 に排他的な CD4-CD8 パラダイムにもかかわらず、αβT 細胞のごく一部 は非従来型 T 細胞で構成されています。これらの非従来型 T 細胞は、 二重陰性 CD4-CD8- T 細胞^{5,12}と二重陽性 CD4+CD8+ T 細胞¹³で構成 されています。非従来型 T 細胞は、最も研究が進んでいない代表的な T 細胞サブセットであり、細胞傷害機能(慢性ウイルス感染)または制御 作用(悪性腫瘍)のいずれかを示すさまざまな状態において重要な役 割を果たすことが確認されています。従来のT細胞と同様に、非従来型

T 細胞も活性化マーカーを発現し、メモリー T 細胞のような特徴を示します¹³⁻¹⁵。CD3+ T 細胞は制御性 T 細胞(Treg)からも構成されており、健康な PBMC 中の CD4+ T 細胞の 3 ~ 5 % を占めます¹⁶。Treg は、CD3、CD4、CD25、CD127 マーカーの相対的発現など、さまざまなアプローチを使って同定されています。Treg 細胞は、胸腺で生成されるナチュラル Treg (nTreg)と末梢誘導性 Treg (iTreg)に細分化されます^{17,18}。Treg 細胞は多くの活性化および記憶特性を示し、自己抗原に対する免疫反応を抑制し、末梢性免疫寛容維持に作用することが知られています^{7,18}。末梢 CD3+ T 細胞の別のサブセットは CD56+ NKT 細胞(リンパ球の 0.04 ~ 1.3 %)であり、自然免疫応答と適応免疫応答の状態に影響を与えます¹⁹。CD16 および CD56 マーカーの発現レベルに基づいて、NKT 細胞はさらにサブセットに類別され、T 細胞と同様に、副刺激分子の関与が NKT 細胞の活性化と働きに影響を与えます^{19,20}。

その他の非従来型 T 細胞には、自然免疫と適応免疫の両方に似た特性 を持ち、自然免疫と適応免疫の橋渡しとなる $\gamma\delta T$ リンパ球があります²¹。 これらの能力により、 $\gamma\delta T$ 細胞は、そのユニークな生物学的特性により、 養子細胞免疫療法の魅力的な候補物質となります。特定の因子とタンパ ク質の発現レベルに基づいて、 $\gamma\delta T$ 細胞はさまざまなサブタイプに細分 化されます²²。

末梢リンパ球には、B 細胞、NK 細胞、ILC などの非 T 細胞も含まれま す^{3,23}。B 細胞は抗体を分泌する CD19+ 発現細胞です。そのため、こ れらの細胞は液性免疫の中心的な要素であり、適応免疫システムの一 部としてほぼすべての各種病原体から保護します。B 細胞は PBMC の 約 15 % を占め、成熟して形質芽球、形質細胞、移行性 B 細胞、メモ リー B 細胞、濾胞性 B 細胞、辺縁帯 B 細胞、制御性 B 細胞、および B-1 細胞に分化することができます。各 B 細胞サブセットは、CD27、 IgM、IgG、IgD、CD38 などのタンパク質と因子のセットの発現に基 づいて特徴付けられます²⁴。PBMC の構成には、病原体に感染した 細胞を直接破壊することで自然免疫システムの重要な構成要素とな る細胞傷害性ナチュラルキラー細胞である NK 細胞(約10%)も含 まれます。CD16 と CD56 マーカーの相対的発現に基づいて、NK 細 胞はさらに異なるサブセットに類別することができます^{3,20}。PBMC は また、系統マーカーと抗原特異的受容体の発現を欠く ILC で構成さ れます。これらの特性にもかかわらず、ILC はサイトカインと分泌タン パク質を生成し、感染した場合の免疫反応の改善につながります。 一部の ILC には、CD4+ T 細胞サブセットと多くの類似点があることが示 されています。したがって、ILC は特定のマーカー発現とサイトカイン産 生に基づいてさらに特定のサブセットに類別できます^{25,26}。

PBMC には DC 細胞(1~2%)も含まれており、これは、ナイーブ CD4+および CD8+T 細胞のプライミングと分化をヘルパーT 細胞、細 胞傷害性エフェクターT 細胞、およびメモリーT 細胞に誘導するとい う独自の特性を持つ専門的な抗原提示細胞として働きます。DC はさ らに、そのオントロジー、局在、表面マーカー発現、サイトカイン産生、 抗原処理、および提示能力に基づいて、骨髄樹状細胞または従来の樹 状細胞(CD1c/cDC2 および CD141/cDC1)と形質細胞様樹状細胞 (CD123/pDC)に類別されます^{27,28}。PBMC にはまた、単球(10~20 %)も含まれ、これは末梢で著しい異質性を示す非常に可塑性の高い自 然免疫細胞タイプです。表面マーカーの相対発現における明確なパターン (CD16、CD14)が、非古典的・中間的・古典的単球という3つの単球 サブセットを区別する基準として認められるようになりました^{3,29}。PBMC はまた、PBMC 分離法に基づいて、好塩基球と呼ばれる細胞の小さな 分画(<1%)でも構成されています。これらの好塩基球は、HLA-DR と CD123 マーカーの相対的発現に基づいて特徴付けられます³⁰。

血液中の各免疫細胞サブセットの複雑な作用を解明することにより、そ れらがどのように免疫反応を形成し、がんや病気の進行、および薬の効 能に影響を与えるかを評価することへの関心が高まっています。この課 題を解決するために、単一細胞レベルおよびサブセットレベルで免疫シス テムのハイスループットかつ詳細な分析を行える方法論と機器が強く求 められています。アジレントは、健康な PBMC における主要な自然免疫 細胞と適応免疫細胞の種類で発現する末梢免疫サブセットの頻度、タン パク質発現、活性化マーカー、疲弊マーカー、および分化マーカーを定 量化するために、45 色スペクトルフローサイトメトリー免疫表現型パネ ルを開発しました。各集団で発現される共通の表面マーカーを使用して 集団頻度を決定し、細胞活性化および増殖(CD69、HLA-DR、CD38)、 細胞疲弊および老化 (PD1、CD223、CD57)、細胞分化 (CCR7、 CD27、CD28、CD45RA、CD45RO、CD127)、細胞可塑性または泳動 電位 (CXCR3、CCR6、CCR5、CXCR5) に関連するマーカーを使用し て各集団のさらなる特性解析を行いました。この詳細な分析は、Agilent NovoCyte Opteon スペクトルフローサイトメータを使用することで可能 になりました。

従来のフローサイトメトリーでは得られる蛍光体のピーク発光が検出器 の数に制限されるが、スペクトルサイトメトリーでは多数の検出器を用い て、すべてのレーザーラインにわたるすべての蛍光体の完全な蛍光スペク トルを捕捉します。そうすることで、スペクトルサイトメトリーでは、一回 の実験で蛍光色素ごとにより多くのパラメータを用いることが可能になり ます^{1,31,32}。

実験方法

試料調製

表 1 は、45 色サンプルを調製するために使用された試薬についての説 明です。

表1. 使用した試薬

試薬	部品番号	メーカー
True-Stain Monocyte Blocker	426102	BioLegend
Brilliant Stain Buffer Plus	566385	BD
CellBlox Blocking Buffer	B001T02F01	Thermo Fisher Scientific
Phosphate Buffer Saline (PBS)	GNM-14190	Genom Biomedical Technology
Fetal Bovine Serum (FBS)	10091148	Thermo Fisher Scientific

染色手順

サンプル前処理法は次のプロトコルに従います。

PBMC の調製

- a. RPMI 1640 完全培養培地を 37 ℃ で少なくとも 30 分間予熱し ます。
- b. PBMC クライオチューブ (約 15 x 10⁶ 個の細胞)を 37 ℃ の ウォーターバスで小さな氷片だけが残るまで解凍します。
- c. クライオチューブの内容物を 50 mL コニカルチューブに移します。
- d. 予熱した 1 mL の RPMI 1640 完全培養培地をクライオチューブに 追加します。ステップ g で使用します。
- e. 50 mL チューブ内の細胞に 5 mL の RPMI 1640 完全培養培地を 一滴ずつ加えます。添加する際は、50 mL チューブを静かに混ぜま す(片手にピペットを持ち、もう一方の手に 50 mL チューブを持 ち、RPMI 1640 完全培養培地を添加しながらチューブを静かに撹 拌します)。
- f. 最初のバッチの 5 mL RPMI 1640 完全培養培地を加えた後は、次の5 mL を少し速く(毎回数滴ずつ)加えます。
- g. 10 mL を加えたら、クライオチューブの内容物を 50 mL チューブに 移します。

- h. 50 mL チューブに RPMI 1640 完全培養培地を追加して 20 mL に します。
- i. 400gで8分間、遠心分離します。
- j. ペレットが動かないように注意して、上澄みを慎重に注ぎ出します。
- k. ペレットを、予熱した RPMI 1640 完全培養培地 2 mL に静かに懸 濁し、20 mL にします。
- I. ステップ i と j を繰り返します。
- m. 後で使用するために、細胞沈殿物を一定量の RPMI 1640 完全培 養培地に懸濁します。

生死判定色素と抗体の混合物の調製

- a. Zombie NIR の生死判定色素を解凍し、0.25 µL をより大きなエッ ペンドルフチューブに移し、1000 µL の 1x PBS (最終希釈 1:4000) を加えます。使用するまで光に当てないようにします。
- b. 4本の1.5 mL エッペンドルフチューブに抗体混合物A、B、C、Dとマークします。優れた染色を行うには、細胞を4段階で染色する必要があります。
- c. 各抗体混合チューブに 10 µL の Brilliant Stain Buffer と 5 µL の True-Stain Monocyte Blocker を加えます。抗体混合チューブ A にのみ 5 µL の CellBlox Blocking Buffer を加えます。
- d. 抗体リスト(表 2)に従って抗体混合物を調製し、A とラベルした チューブにグループ A の抗体を加え、B とラベルしたチューブにグ ループ B の抗体を加える、というように繰り返します。
- e. 混合物チューブを16,000 g ~ 18,000 g で 5 分間遠心分離します。 抗体を使用する場合は、液体の上部から混合物を吸引し、チューブ の底に触れないようにしてください。抗体の凝集体が底に残ります。

表 2. 抗体染色グループ (CD45RO SBUV445 抗体を除く)

	特異性	蛍光色素	添加量(µL)
染色	グループ A		
1	CCR6	BV711	2.5
2	CCR5	BUV563	2.5
3	TCRγδ	PerCP-eFluor 710	0.63
4	CXCR5	BV750	0.63
5	IgG	BV605	0.63
6	CXCR3	PE-Cy7	1.25
7	CD223 (LAG-3)	NovaFluor Blue 660/120S	1.25
8	CD69	StarBright Blue 765	1.25
9	CCR7	BV421	2.5
染色	グループ B		
1	CD20	Spark YG 593	1.25
2	CD1c	Alexa Fluor 647	2.5
3	CD28	BV650	5
4	PD-1	BV785	5
5	CD159c	PE	10
染色	グループ C		
1	CD127	Spark Red 718	1.25
2	CD2	PerCP-Cy5.5	2.5
3	CD337	PE-Dazzle 594	2.5
4	CD3	BV510	2.5
5	CD27	APC-H7	2.5
6	CD25	PE-Alexa Fluor 700	5
7	CD11c	eFluor 450	5
染色	グループ D		
1	CD45RA	BUV395	0.32
2	IgD	BV480	0.32
3	CD4	CF594	0.32
4	CD14	Spark Blue 550	0.32
5	HLA-DR	PE-Fire 810	0.32
6	CD33	StarBright Blue 580	0.32
7	CD16	BUV496	0.63
8	CD57	FITC	0.63
9	CD24	PE-Alexa Fluor 610	0.63
10	CD95	PE-Cy5	0.63
17	CD38	APC-Fire 810	0.63
12	CD31		0.63
13	CD141	BB315	1.25
14	CD8		1.20
16	CD159a	APC.	2
17	CD314	BUV615	2.5
18	CD56	BUV737	2.5
19	CD45	PerCP	2.5
20	IgM	BV570	2.5
21	CD39	BUV661	2.5
22	CD123	Super Bright 436	2.5
			·J

染色

- a. FACS チューブをマークします。未染色の対照群チューブ、単一染色 チューブ、およびマルチカラーチューブをマークします。
- b. 各マルチカラーチューブに細胞懸濁液 100 μL(約3 x 10⁶ 個の細胞)を加え、各未染色の対照群チューブと単一染色チューブに細胞 懸濁液 10 μL(約3 x 10⁵ 個の細胞)を加えます。
- c. マルチカラーチューブと生死判定用単一染色チューブを PBS で洗 浄し、その他の単一染色チューブと未染色の対照群チューブを洗浄 バッファ (PBS + 2 % FBS) でチューブあたり 3 mL で洗浄します。
- d. 500gで5分間遠心分離し、上澄みを慎重に注ぎ出します。
- e. 100 μL の Zombie NIR 希釈液をマルチカラーチューブと生死判定 用単一染色チューブに加え、よくボルテックスし、暗所で 30 分間氷 上に置きます。
- f. 同時に、他の単一染色チューブを表 2 に示す最適な抗体レベルで 染色し、暗所で 30 分間氷上に置きます。anti- CD223 NovaFluor Blue 660/120S シングル染色チューブに CellBlox Blocking Buffer を追加します。
- g. 30 分間インキュベーションした後、マルチカラーチューブ、生死判 定用単一染色チューブ、およびその他の単一染色チューブを 3 mL の洗浄バッファで洗浄します。
- h. ステップ d を繰り返します。
- i. 洗浄バッファ 200 µL を生死判定用単一染色チューブ、未染色の対 照群チューブ、およびその他の単一染色チューブに追加し、ボルテッ クスで混合して氷上で保存します。
- j. すべてのマルチカラーチューブに anti-CD45RO SBUV445 を加え、 ボルテックスし、暗所で 30 分間氷上に置きます。
- k. 洗浄せずに、抗体混合物 A をすべてのマルチカラーチューブに加え てボルテックスし、暗所で30 分間氷上に置き、その後室温(RT)で 20 分間置きます(これにより、CCR および CXCR マーカーの染色 が改善されます)。
- すべてのマルチカラーチューブを洗浄バッファ 3 mL で 1 回洗浄します。
- m. ステップ d を繰り返します。

- n. すべてのマルチカラーチューブに抗体を加えてボルテックスし、暗所 で 40 分間氷上に置きます。
- すべてのマルチカラーチューブを洗浄バッファ 3 mL で 1 回洗浄し ます。
- p. ステップ d を繰り返します。
- q. マルチカラーチューブに抗体混合物 C を加えてボルテックスし、暗 所で 40 分間氷上に置きます。
- r. すべてのマルチカラーチューブを洗浄バッファ 3 mL で 1 回洗浄します。
- s. ステップ d を繰り返します。
- t. マルチカラーチューブに抗体混合物 D を加えてボルテックスし、暗 所で 40 分間氷上に置きます。
- u. すべてのマルチカラーチューブを洗浄バッファ 3 mL で 2 回洗浄し ます。
- v. ステップ d を繰り返します。
- W. 洗浄バッファ 300 µL をマルチカラーチューブに加えてボルテックスし、氷上で保存します。

結果と考察

抗体価測定

このアプリケーションノートでは、この 45 色パネル(表 3 に概要を示し ます)で提案されているすべてのマーカーと蛍光色素の組み合わせが、 抗体価測定分析によって評価されています。図 1 に示す各抗体の必要量 は、次のように計算される Stain index から決定されます。(MFI (陽性) -MFI (陰性)) / (2xSD (陰性))、ここで MFI は平均蛍光強度です。最適 な抗体濃度は、Stain index が最大または最大に近い濃度か、あるいは 拡散誤差がより低くなるような、最低濃度が選択されます。

このパネル設計の主な目的は、以前に公開された 40 マーカーパネル である OMIP-069 を元に、さらに高性能なものを構築することでした。 OMIP-069 パネルに加えて、さらに CD45RO、CD33、CD233 (LAG-3)、 CD69、CD31 の 5 つのマーカーも含めました。抗体と蛍光色素の全パ ネルを表 3 に示します。立体障害によって誘発されるこのような大きな パネルでの染色は複雑な性質となるため、抗体-抗原結合を最適化して 4 つの染色グループを特定し、サンプルとともに連続パターンでインキュ ベートしました (表 2)。

表 3. 抗体情報

	マーカー	蛍光色素	クローン	メーカー	部品番号
1	CD45RA	BUV395	5H9	BD Biosciences	740315
2	CD45RO	StarBright UltraViolet 445	UCHL1	Bio-RAD	MCA461SBUV445
3	CD16	BUV496	3G8	BD Biosciences	612945
4	CCR5 (CD195)	BUV563	2D7/CCR5	BD Biosciences	741401
5	CD314 (NKG2D)	BUV615	1D11	BD Biosciences	751232
6	CD39	BUV661	TU66	BD Biosciences	749967
7	CD56	BUV737	NCAM16.2	BD Biosciences	612766
8	CD8	BUV805	SK1	BD Biosciences	612889
9	CCR7 (CD197)	BV421	G043H7	BioLegend	353208
10	CD123	Super Bright 436	6H6	Thermo Fisher Scientific	62-1239-42
11	CD11c	eFluor 450	3.9	Thermo Fisher Scientific	48-0116-42
12	IgD	BV480	IA6-2	BD Biosciences	566187
13	CD3	BV510	SK7	BioLegend	344827
14	lgM	BV570	MHM-88	BioLegend	314517
15	lgG	BV605	G18-145	BD Biosciences	563246
16	CD28	BV650	CD28.2	BioLegend	302945
17	CCR6 (CD196)	BV711	G034E3	BioLegend	353435
18	CXCR5 (CD185)	BV750	RF8B2	BD Biosciences	747111
19	PD-1 (CD279)	BV785	EH12.2H7	BioLegend	329929
20	CD141	BB515	1A4	BD Biosciences	565084
21	CD57	FITC	NK-1	BD Biosciences	555619
22	CD14	Spark Blue 550	63D3	BioLeaend	367148
23	CD33	StarBright Blue 580	WM53	Bio-RAD	MCA1271SBB580
24	CD223 (LAG-3)	NovaFluor Blue 660/120S	3DS223H	Thermo Fisher Scientific	H048T02B08
25	CD45	PerCP	HI30	Agilent Technologies	8931017
26	CD2	PerCP-Cv5.5	TS1/8	BioLeaend	309225
27	ΤCRγδ	PerCP-eFluor 710	B1.1	Thermo Fisher Scientific	46-9959-41
28	CD69	StarBright Blue 765	FN50	Bio-RAD	MCA2806SBB765
29	CD31	StarBright Blue 810	WM59	Bio-RAD	MCA1738SBB810
30	CD159c (NKG2C)	PE	134591	R&D Systems	FAB138P-100
31	CD20	Spark YG 593	2H7	BioLegend	302367
32	CD337 (Nkp30)	PE-Dazzle 594	P30-15	BioLegend	325231
33	CD4	CF594	C4/206	Biotium	#BNC940206-500
34	CD24	PE-Alexa Fluor 610	SN3	Thermo Fisher Scientific	MHCD2422
35	CD95 (FAS)	PE-Cy5	DX2	BioLegend	305610
36	CD25	PE-Alexa Fluor 700	CD25-3G10	I hermo Fisher Scientific	MHCD2524
20		DE Eiro 910	GUZ3H7	Piel egond	207692
39	CD159a (NKG2A)	APC	RFA110	Miltenvi	130-113-563
40	CD1c	Alexa Fluor 647	L161	BioLegend	331510
41	CD19	Spark NIR 685	HIB19	BioLegend	302269
42	CD127	Spark Red 718	A019D5	BioLegend	351376
43	Viability	Zombie NIR	N/A	BioLegend	423105
44	CD27	APC-H7	M-T271	BD Biosciences	560223
45	CD38	APC-Fire 810	HB-7	BioLegend	356643

A. UV レーザー (349 nm)

D. 黄レーザー (561 nm)



図 1.45 色パネルの抗体価測定。すべての抗体は、健康な個人から採取した白血球パックから分離し凍結保存された PBMC を使用して、染色量 100 μL あたり 1 ~ 20 μL の 範囲の開始濃度で 6 段階の濃度で滴定されます。X 軸は希釈系列を示し、6 つの希釈サンプルはすべて同じプロットに表示されています。赤い四角は、パネルに使用される最 終的に選択された抗体濃度を強調表示したものです。最適な抗体濃度は Stain index (Stain index が最大または最大に近い濃度か、あるいは拡散誤差がより低くなるような 最低濃度) に基づいて選択され、赤いボックスで示されています。測定データはすべて、Agilent NovoExpress 2.0.0 ソフトウェアによって分離されました。デブリとダブレットは 除外され、ファイルは FlowJo v10.10.0 を使用して連結されました。(A) 349 nm レーザー励起蛍光色素、(B) 405 nm レーザー励起蛍光色素、(C) 488 nm レーザー励起 蛍光色素、(D) 561 nm レーザー励起蛍光色素、(E) 637 nm レーザー励起蛍光色素。Stain index の計算は次のとおりです。(MFI (陽性) -MFI (陰性)) / (2xSD (陰性))、 ここで MFI は平均蛍光強度です。

図 2 に示されているように、自己蛍光 (AF) 減算用の 2 つの集団と共 に、使用したマーカーごとにリファレンススペクトルコントロールが実行さ れました。

A. 45 種類の蛍光色素のスペクトルシグネチャ





図 2.45 色のパネルデータをアンミックス するために使用されたスペクトルシグネ チャ (45 + 2 AF)。単一染色の対照群は、 Agilent NovoCyte Opteon スペクトルフ ローサイトメータを用いてデフォルトのゲイン 設定で取得されたものです。蛍光色素のスペ クトルシグネチャ (正規化された発光スペク トル)は、一次レーザーと自己蛍光によって グループ化されたスペクトルプロットとして 表示されています。(A) 45 個の蛍光色素、 (B) 2 つの自己蛍光シグネチャ パネルをさらに最適化するために、Agilent NovoExpress ソフトウェアを 使用して、類似性指数と Spillover Spreading マトリックスを計算しました (図 3 および 4)。

CD4ERA RUV20E	1																																												
CD45RA BUV395	0.70	1																																											
CD16 BUV/96	0.73	0.47	1																																										
CCP5(CD195) BLIV563	0.23	0.17	0.2	1																																									
CD314(NKG2D) BLIV615	0.15	0.08	0.12	0.51	1																																								
CD39 BLIV661	0.00	0.00	0.02	0.00	0.27	1																																							
CD56 BUV737	0.05	0.05	0.03	0.03	0.37	0.20	1																																						
CD8 BUV805	0.05	0.00	0.05	0.05	0.05	0.00	1 26	1																																					
CCP7(CD197) BV/421	0.33	0.32	0.05	0.00	0.03	0.03	0.33	0.05	1																																				
CD123 Super Bright 436	0.11	0.25	0.03	0.02	0.01	0.01	0.01	0.05	1 92	1																																			
CD11c eFluor 450	0.09	0.32	0.15	0.05	0.03	0.01	0.01	0.04	0.74	0.02	1																																		
InD BV/80	0.03	0.52	0.13	0.05	0.04	0.01	0.01	0.04	0.23	0.38	151	1																																	
CD3 BV510	0.02	0.10	0.54	0.00	0.16	0.05	0.01	0.02	0.12	0.22	0.2	0.8	1																																
IgM BV570	0.03	0.06	0.00	0.49	0.10	0.03	0.04	0.04	0.12	0.20	1.25	0.2	142	1																															
laG BV605	0.03	0.07	0.00	0.32	0.54	0.23	0.08	0.04	0.1	0.11	01 0	112	138 0	73	1																														
CD28 BV650	0.03	0.07	0.04	0.06	0.24	0.41	0.00	0.04	0.11	0.11	0.1	0.12	0.2 0	21 0	162	1																													
CCP6(CD196) BV711	0.03	0.05	0.04	0.00	0.08	0.28	0.10	0.11	0.02	0.02	0.03	2.00	112 0	14	0.2 (150	1																												
CXCR5(CD185) BV750	0.03	0.05	0.04	0.02	0.00	0.07	0.4	0.14	0.00	0.00	103	0.03	106 0	06 0	116 0	127	0.6	1																											
PD_1(CD279) BV785	0.02	0.01	0.01	0.01	0.03	0.04	0.23	0.23	0.04	0.08	0.03	103		06 0	11 (17	0.0	0.77	1																										
CD141 BB515	0.02	0.04	0.14	0.11	0.03	0.04	0.01	0.01	0.00	0.01	0.07	0.03	103 0	02 0	0.01	0	0.00	0	0	1																									
CD57 FITC	0.02	0.01	0.13	0.14	0.02	0	0.01	0.01	0	0.01	0.01	0.08	1 05 0	04	0.02	0	0	0	ŏ	0.98	1																								
CD14 Spark Blue 550	0.01	0.01	0.13	0.14	0.02	0.01	0.01	0.01	0.01	0.01	102	0.00	113 0	16	0.1 0	04	0.03	0.02	0.01	0.50	0.77	1																							
CD33 SBB580	0.01	0.01	0.11	0.55	0.24	0.05	0.00	0.01	0.01	0.01	1.02	01	1 26 0	58	0.4 0	112	30.0	0.02	0.02	0.22	0.31	0.59	1	1																					
CD223(LAG-3) NEBlue 660/120s	0	0.01	0.01	0.05	0.05	0.43	0.24	0.03	0	0.01	0 0	2.01	0.01 0	05 0	0.4 0	117	0.17	0.03	0.02	0.00	0.01	0.55	0.17	1																					
CD45 PerCP	0	0.01	0.02	0.07	0.28	0.45	0.24	0.04	0	0.01 (0.01	0.01	108 0	16 0	34 0	152	0.33	0.19	0.02	0.00	0.02	0.08	0.13	0.53	1																				
CD2 PerCP-Cv5 5	0	0.01	0.01	0.05	0.22	0.44	0.42	0.08	0	0.01	0.01	0.01	0.08 0	15 (32 0) 51	0.61	0.3	0.19	0.01	0.02	0.09	0.13	0.57	0.76	1																			
TCRv& PerCP_eEluor 710	0.01	0.01	0.02	0.03	0.14	0.33	0.5	0.09	0	0.01	0.01	0.01	106	1 (123 (38	0.65	0.32	0.2	0.01	0.02	0.08	0.11	0.46	0.47	0.9	1																		
CD69 SBB765	0.02	0.01	0.02	0.06	0.06	0.09	0.5	0.05	0.01	0.02	0.01	0.01	0.00	08 0	14 () 17	0.39	0.52	0.54	0.01	0.02	0.00	0.11	0.40	0.26	0.46	0.52	1																	
CD31 SBB810	0.02	0.04	0.00	0.00	0.04	0.05	0.40	0.28	0.01	0.02	0.02	0.00		06 0	000 0	11	0.24	0.33	0.54	0.00	0.07	0.14	0.14	0.14	0.15	0.40	0.34	0.83	1																
CD159c(NKG2C) PE	0.00	0.01	0.02	0.61	0.28	0.05	0.01	0.01	0	0 0	0.01	0.03	1.08 0	64 (32 (0.05	0.02	0.1	0.00	0.07	0.11	0.26	0.67	0.08	0.08	0.07	0.05	0.05	0.05	1															
CD20 Spark YG 593	0	0	0.01	0.46	0.61	0.22	0.04	0.01	0	0	0	0	0 0	47 (142 (112	0.02	0	0	0.01	0.02	0.07	0.27	0.00	0.24	0.25	0.19	0.06	0.04	±	1														
CD337(NKp30) PE-Dazzle 594	0	0	0.02	0.37	0.64	0.18	0.05	0.01	0	0 (011	011	105 0	12 (148 (16	0.05	0.02	0.01	0.04	0.06	0.19	0.43	0.24	0.31	0.20	0.21	0.00	0.08	0.52	0.70	1													
CD4 CE594	0	0	0.02	0.27	0.71	0.22	0.03	0.01	0	0	0	0 0		33 (143 (115	0.03	0.01	0.01	0.04	0.00	0.03	0.15	0.07	0.26	0.23	0.16	0.03	0.02	0.35	0.83	88.0	1												
CD24 PE-Alexa Eluor 610	0	0	0.01	0.21	0.57	0.33	0.00	0.01	0	0 (011	201	103 0	20 0	141 0	127	0.13	0.04	0.02	0.02	0.04	0.13	0.26	0.42	0.20	0.55	0.44	0.16	0.12	0.00	0.60	0.84	0.78	1											
CD95 PE-Cv5	0.01	0.01	0.01	0.09	0.37	0.46	0.18	0.03	0	0 0	0.01	0.01	0.02 0	15 (28 (28	0.15	0.05	0.03	0.01	0.02	0.08	0.14	0.59	0.68	0.68	0.52	0.2	0.13	0.14	0.53	0.49	0.5	183	1										
CD25 PE-Alexa Eluor 700	0	0	0	0.09	0.22	0.23	0.31	0.04	0	0	0	0 1	0.01 0	13 (18 (15	0.21	0.11	0.05	0.01	0.02	0.07	0.12	0.26	0.35	0.64	0.74	0.38	0.23	0.15	0.41	0.29	0.3	152	0.59	1									
CXCR3(CD183) PE-CV7	0.01	0.01	0.01	0.02	0.06	0.06	0.16	0.11	0	0	0	0 1	0.01 0	03 0	0.05 (0.04	0.07	0.1	0.12	0	0	0.02	0.03	0.06	0.1	0.22	0.25	0.52	0.52	0.02	0.13	0.07	0.08	0.14	0.18	137	1								
HLA-DR PE-Fire 810	0	0	0	0.12	0.1	0.06	0.11	0.1	0	0	0	0 0	0.01 0	12 (0.09 0	0.04	0.05	0.06	0.09	0.01	0.02	0.06	0.13	0.07	0.1	0.2	0.23	0.37	0.54	0.2	0.24	0.16	0 1 4	19	0.18	0.31	0.74	1							
CD159a(NKG2A) APC	0	0	0.01	0.04	0.21	0.8	0.3	0.06	0	0	0 0	0.01	0.02 0	09 (18 (0.36	0.26	0.07	0.04	0	0	0.01	0.03	0.54	0.34	0.39	0.34	0.08	0.04	0.06	0.26	0.2	0.26	0.36	0.54 0	26	0.08	0.07	1						
CD1c Alexa Fluor 647	0	0	0	0.01	0.07	0.72	0.3	0.05	0	0	0	0	0 0	02 (0.04	122	0.25	0.03	0.02	0	0	0	0.01	0.56	0.25	0.34	0.3	0.05	0.02	0.03	0.1	0.07	0.1	0.16	0.34	14	0.03	0.03	0.88	1					
CD19 Spark NIR 685	0	0	0	0.01	0.06	0.63	0.34	0.07	0	0	0	0	0 0	02 (0.05 (21	0.31	0.05	0.03	0	0	0	0.01	0.47	0.25	0.4	0.36	0.07	0.03	0.02	0.08	0.06	0.09	0.14	0.31	0.17	0.05	0.04	0.72	0.92	1				
CD127 Spark Red 718	0.08	0.11	0.09	0.04	0.04	0.45	0.57	0.19	0.02	0.03	0.03	0.02	0.05 0	02 0	0.04 () 15	0.4	0.1	0.07	0.01	0	0.01	0.01	0.3	0.12	0.33	0.42	0.13	0.06	0.01	0.03	0.02	0.03	0.07	0 17 0	0.18	0.06	0.04	0.52	0.62	0.73	1			
L/D Zombie NIR	0	0	0	0	0.02	0.23	0.6	0.22	0	0	0	0 1	0.01 0	.01 (0.03	0.08	0.27	0.21	0.14	0	0	0.01	0.01	0.2	0.12	0.26	0.33	0.34	0.18	0.01	0.04	0.03	0.03	0.07	0.15	0.23	0.21	0.1	0.33	0.35	0.41	0.67	1		
CD27 APC-H7	0	0	0	0	0.02	0.1	0.31	0.42	0	0	0	0	0 0	.01 (0.02 (0.04	0.14	0.14	0.19	0	0	0	0	0.07	0.04	0.13	0.16	0.23	0.18	0.01	0.03	0.01	0.02	0.04	0.07	0.11	0.3	0.23	0.15	0.16	0.21	0.41	0.64	1	
CD38 APC-Fire 810	0	0	0	0	0.02	0.11	0.24	0.44	0	0	0	0	0 0	.01 (0.02	0.04	0.11	0.09	0.16	0	0	0	0	0.07	0.04	0.1	0.13	0.18	0.19	0.01	0.03	0.02	0.02	0.04	0.07	0.08	0.2	0.27	0.16	0.15	0.18	0.34 0	0.38	0.8	1
Complexity Index: 55.38	45RA BUV395	45RO SBUV445	16 BUV496	35(CD195) BUV563	314(NKG2D) BUV615	39 BUV661	56 BUV737	8 BUV805	37(CD197) BV421	123 Super Bright 436	11c eFluor 450	BV480	3 BV510	1 BV570	BV605	28 BV650	36(CD196) BV711	CR5(CD185) BV750	-1(CD279) BV785	141 BB515	57 FITC	14 Spark Blue 550	33 SBB580	223(LAG-3) NFBlue 660/120s	45 PerCP	2 PerCP-Cy5.5	ኛላδ PerCP-eFluor 710	59 SBB765	31 SBB810	159c(NKG2C) PE	20 Spark YG 593	337(NKp30) PE-Dazzle 594	4 CF594	24 PE-Alexa Fluor 610	95 PE-Cy5	25 PE-Alexa Fluor 700	CR3(CD183) PE-Cy7	A-DR PE-Fire 810	159a(NKG2A) APC	1c Alexa Fluor 647	19 Spark NIR 685	127 Spark Red 718	Zombie NIR	27 APC-H7	38 APC-Fire 810
	8	9	9	0 S	8	8	8	0	S	9	8	ЪВ	8	2B	96	8	S	Š	9	8	8	8	8	9	9	9	1C	8	0	8	8	8	8	8	0	8	Š	Ŧ	8	8	8	8	5	8	8

図3.45 色のパネルで使用される蛍光色素の類似性指数マトリックス

Source\Target	CD45RA BUV395	CD45RO SBUV445	CD16 BUV496	CCR5(CD195) BUV563	CD314(NKG2D) BUV615	CD39 BUV661	CD56 BUV737	CD8 BUV805	CCR7(CD197) BV421	CD123 Super Bright 436	CD11c eFluor 450	IgD BV480	CD3 BV510	IgM BV570	IGG BV605	CD28 BV650	CCR6(CD196) BV711	CXCR5(CD185) BV750	PD-1(CD279) BV785	CD141 BB515	CD57 FITC	CD14 Spark Blue 550	CD33 SBB580	CD223(LAG-3) NFBlue 660/120s	CD45 PerCP	CD2 PerCP-Cy5.5	TCRyð PerCP-efluor 710	CD69 SBB765	CD31 SBB810	CD159c(NKG2C) PE	CD20 Spark YG 593	CD337(NKp30) PE-Dazzle 594	CD4 CF594	CD24 PE-Alexa Fluor 610	CD95 PE-Cy5	CD25 PE-Alexa Fluor 700	CXCR3(CD183) PE-Cy7	HLA-DR PE-Fire 810	CD159a(NKG2A) APC	CD1c Alexa Fluor 647	CD19 Spark NIR 685	CD127 Spark Red 718	L/D Zombie NIR	CD27 APC-H7	CD38 APC-Fire 810
CD45RA BUV395		11.1	8.2	2	0.9	0.8	1	0.9	1.8	2.8	2.7	1	2.5	0.7	0.3	0.7	0.8	0.7	0.5	1.7	1.6	1.2	1.6	0.4	1	2.3	1.1	0.3	0.5	2.3	0.8	0.8	1.2	0.5	0.3	0.9	0.4	0.8	0.7	0.7	0.7	0	0.3	0	0.7
CD45RO SBUV445	2.5		9	2.4	1.2	0.8	1.1	1	2.9	4.5	3.9	2.3	2.2	0.7	0.4	0.8	0.6	0.5	0.4	1.7	2.9	0.8	1.8	0.3	1.1	1.9	1.2	0.1	0.6	2.3	0.8	0	1.7	0.5	0.6	0.6	0.5	0.8	1	0.8	1.1	0.4	0.3	0	0.6
CD16 BUV496	1.5	81		2	0.9	0.6	1	1	11	0.9	1	11	24	1.5	0	0.7	0.5	12	0	51	5.5	29	13	0.7	0.8	1.5	1	0	0	22	19	1.8	21	16	19	0.7	0	0.6	0.5	0.2	17	0	0.9	0	12
CCR5(CD195) BLIV563	1.8	8	53		1.4	1.4	15	0.7	11	0.5	23	1.5	23	1.5	0.9	1.4	17	1.7	0.5	8	0.3	3.5	4.8	1.6	2.8	23	26	11	0	67	1.0	5.3	53	3.0	3.4	1.8	21	0.8	3.1	5.2	1.4	1.6	1	13	0.3
CD214(NKC2D) BUV615	1.0	10	21	2	1.44	2.5	2.6	2.1	0	22	1.4	0.0	2.0	1.0	1.0	2.5	2.2	27	2.1	5.5	9.0	2	4.0	2.6	2.0	5.1	5.9	2.5	2.2	6.5	Q./	6.0	77	6.0	6.5	27	2.1	2.5	5	6.7	4.7	2.1	1.4	0.0	11
CD20 BUV661	1.9	5	4	21	1.0	2.0	2.0	2.1	0	2.4	1.4	1.5	2.0	1.9	0	2.0	2.4	2.7	1.4	2.2	0	21	4.2	2.0	2	5.7	6.7	2.0	2.2	47	6.6	7.6	21	6	0.0	2.0	27	2.0	12.2	20.6	11.1	5.6	4.4	1	2.0
CD56 BUV727	1.2	0	22	1.1	0.7	1.4	0.9	2.4	1.5	1.9	27	1.3	2.2	0.9	0.6	0.0	2.4	2.0	1.4	0.0	0	0.1	2.4	0.9	2.0	5.0	12	6	1	2.2	2.5	27	0.0	2.0	1.5	3.2	2.7	2.5	4.4	10.2	0.0	9.7	10	7.0	7.6
CD9 PLIV905	20	6.0	2.0	1.1	0.7	0.7	16	2.0	1.0	1.0	1	0.6	1.0	0.0	0.0	1.1	0	0.7	9.2	0	0	0.1	1.5	0.0	1.1	2.5	1.0	2.2	27	17	1.0	0.6	1.4	1.6	0.6	1.0	2.2	2.0	0.0	0.2	1.7	1.6	2.1	7.0 1	11.7
CD8 B0 V805	1.1	3.3	0.1	1.1	0	0.7	1.5	0.2	- 4	7.0	E 0	1.0	1.9	0.3	0.6	0	0	2	0.0	0.2	2	1	0	0.7	0	2.0	0	2.3	0.7	1.7	0	0.0	0	0	0.0	0	0	0.4	0.0	2	0	0	2.1	0.6	0.7
CCR7(CD197)BV421	1.0	0.2	0.0	1.1	0	0	0.5	0.5		7.9	3.0	1.9	1.0	0	0.0	0	0	0	0	1.0	0	0.7	1	0	0.7	1.0	0	0	0	1.9	0	0	0	0	0	0	0	0.4	1	1	1.7	0	0	0.0	0
CD125 Super Bright 450	1.0	0.0	2.7	0.0	0.0	0.4	1	0.5	4.4	10.4	1.2	4.4	4.0	1.6	1	1.7	0	0	0.0	1.0	0	2.7	1.4	0.5	0.7	0	1.1	0	0.0	2.2	1.1	0	0.0	1.4	0.4	0.4	0.0	0	1.1	0.4	1.7	1	0	0.4	1
CD11C 01000 450	2.0	2.0	47	1.4	0.2	0.4	0	0.4	0.7	12.4	0.0	4.Z	4.2	1.0	1.0	0.7	0.7	0	0.0	1.9	4.0	2.4	1.4	0.0	0.4	0.1	0	0	0.0	1.7	0	0.0	0.5	1.4	0.4	0.4	0.9	0.2	0	2.4	0	0	0	0.0	0.7
IGD 8V460	1.2	4.8	4.7	1.4	1.6	1.0	1.0	1	2.5	3.8	2.8	0.4	1.9	1.9	1.2	1.0	0.7	0.1	1.7	4.4	4.0	1.4	1.3	0.0	0.1	2.1	0	0.7	0	1.7	1.0	1.7	1.6	1.6	0.1	0.7	0	0.3	1.6	1.0	1.6	0.0	0.7	0	J./
LD3 BV510	1.0	0.9	5.3	2.3	1.0	1.2	1.Z	0	2.3	3.Z	2.5	2.4	4.1	2.5	2.3	1.8	2	2.1	0.0	3	3.5	1.7	1.0	0.3	2.1	2.5	2.3	0.7	0.9	2.0	1.8	4.0	1.0	1.0	2.1	0.7	0	1.0	1.0	1.3	1.0	0.8	0.7	0.3	J./
IGINI BV570	1.2	2.4	7.0	3.Z	1.5	3.9	0.7	0.0	0.8	10	0	1.6	4.1	0	2	0	3.0	0	2.0	0.0	0	10	0.7	1.6	4.9	5.9	6.1	0	2.2	4.4	3.5	4.3	0.7	3.4	0.3	0	2.9	1.3	1.1 A.E.	0	0	2.1	1.0	1.0	-
IGG BV605	0	3.5	7.2	5.1	1.9	1.9	1.0	1.5	0.7	13	0	1.0	2.5	3	0.0	0	2.0	3.3	2.5	0	1.0	1.3	2.7	1.0	4.4	5.Z	0.1	2.0	1.7	3.2	4.1	5.0	0.7	7.3	6.5	2.3	3.0	5.0	4.5	10.0	3.5	4	1.6	0.0	0
CD26 BV650	0.2	3	2.2	0.7	0.6	2.2	1.8	1.5	2.7	2.5	2.8	0.3	17	0.7	0.8	1.0	4.1	5.9	3	0	1.9	0.3	2.3	1.0	4	7.1	7.0	1.9	1.7	3.7	5.Z	0.7	2.0	3.2	0.4	2.8	2	1.8	6.1	12.3	7.9	4	2.3	2.8	2.5
CCR0(CD190) BV / 11	0.8	3.0	2.0	1	0	0.0	2.2	1.5	2.0	2.3	2.4	0	0	0.9	0	1.3	0.0	5.7	4.5	0	0	0	0	0	3.4	7.1	7.5	2.7	2.9	1.9	2.0	3.7	1.9	3.0	0	2.0	2.3	1.5	0.1	12.1	9.4	0.7	3	4.1	3.0
CACR5(CD165) BV / 50	0	1.0	0	1	0	0.8	2	1.7	1.9	0	1.1	0	0	0	0.6	0.7	2.8	C 4	5.8	0	0	0	0	0	2.9	2.5	3	3.4	2.7	2.0	1.9	0	1.2	1.9	0	0.2	1.8	1.5	0	3.4	2.5	3.3	2.9	3.Z ·	3.9
PD-1(CD2/9) BV/05	0	0	0.0	10	0.6	0.0	1.2	1.0	2.3	1.0	1	0.5	0	0	0	0.7	1.0	3.4		1.7	10.0	1.0	10	0.4	1.0	0.0	0.4	2.9	3.0	0	0.5	1.2	0	0.9	0.0	0.7	2.0	2.1	1.1	2.1	1.3	1.9	2	0.0	4.Z
CD141 BB515	1.3	0	2.6	1.Z	0	0.4	0.3	0.3	0.6	0	1.8	1.5	0	0.6	0	0.4	1.6	1.7	0	10.0	19.9	0.1	1.2	0	0.2	2.2	3.4	0.4	0	0	0	07	1.4	0	2.1	0.7	0.8	1.2	1.Z	0	2.1	0.5	0	0	0
CD57 FITC	0.3	2	1.4	0.3	0.3	0.2	0.2	0.2	0.5	0.4	0.0	0.6	1.3	1.3	0.4	0.5	0.7	0.9	0.0	19.9	40.0	8.5	1.5	0.6	0.8	1.0	1.4	0.7	0.5	0.9	0.6	2.7	1.5	1.5	1.2	0.5	0.3	0.4	0.3	1.Z	0.8	0.6	0.2	0.2	J.3
CD14 Spark Blue 550	0	2.5	1.6	0.6	0	0.5	0.3	0.2	1.3	1.2	0.8	0.5	1.6	1.6	0.9	0.8	1.1	1.2	0.9	12.7	19.8	4.0	2.9	0.9	1.3	2.3	3.2	1.2	1.4	70	1.4	3.5	2.4	2.1	2.1	1.2	0.8	0.6	1.8	3.Z	0.7	0.7	0.5	0	J.6
CD33 SBB580	0	2.1	1.4	5.3	1.1	0.8	0.5	0.5	2.7	0	1.7	0.5	0	5.5	1.9	1.1	1.2	1.2	0.9	3.9	5.5	4.8		1.3	2	2.9	3.7	1.4	1.1	7.9	3	5.8	3.4	3.6	3	1.5	1.2	0.8	2.3	3.1	2.2	1.1	0.4	0	0
CD223(LAG-3) NFBlue 660/120s	0	2.1	2	0.5	0.4	0.8	1.1	0.7	1.3	2.2	0.3	0	0.4	1.2	0	2.5	2.8	2.5	1.4	6	6.1	1.9	2.2		3.3	6.4	6.5	3.5	3.2	3.4	4.4	4.8	2.6	5.2	4.2	2.4	2.2	1.4	9.9	16.9	8.9	5	2.6	3.3	3.6
CD45 PerCP	0.3	2.3	2.2	1.1	0.6	2	2.4	1.5	1.5	2.5	2	0.5	1.8	1	0.7	2.1	3.6	4	2.8	2.1	2.7	0.9	3.7	3		6.2	7.3	3.5	3	4.5	6.9	7.5	2.9	7.1	9.1	3.3	2.5	2.4	6.4	10.7	7.8	3.5	2.2	2.5	2.6
CD2 PerCP-Cy5.5	0	0	2.9	0.8	0	1.6	2.4	2	2	2.2	2.5	0.3	1.2	0.9	1.1	2.4	6.4	6	4.5	3.2	2.5	0	3.4	2.5	6.5		12.1	5.6	4.4	4.5	6	5.9	2.6	7.1	/	4.8	4	3.2	7.8	15.2	12	6.1	3.1	4	4
TCRyo PerCP-eFluor 710	1.2	5.2	0	0	0.4	0.7	1.9	1.1	1.9	3.2	0	0	1.2	0.9	0.3	1.3	7.6	6.8	6.6	3.6	6.7	0	0	2.3	5.2	16		3.1	0	3.6	4./	0.6	0	4.4	4.2	5.7	2.6	3.1	5.9	7.1	10	5.9	3.2	0	3.8
CD69 SBB/65	0	0	6.9	1.2	0	1.2	4.3	1.1	2.9	14.7	4.3	0	1.3	0	1.9	0	1.5	5.8	8.2	0	0	1.8	0	1.3	0	0	3.5		4./	4.5	0	5.5	0	0	4.4	2.7	1.7	3.8	0	0	4.5	0.5	4.6	5.1	5.2
CD31 SBB810	1.5	3	1.8	1.4	1.1	1.2	0.8	2.6	2.2	0.9	1.6	0.5	4.6	0.9	1.1	0.8	3.2	6./	6.8	4.8	9.1	7.4	5.4	0.8	2.6	6.2	4.1	17		1.5	2.1	4	2.8	1.6	1.1	1.5	1.7	2.6	3	8.9	2.2	3.1	1.4	3.2	0.4
CD159c(NKG2C) PE	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0		0	0	0	0	0	0	0	0	0	0	0	0	0	0	0
CD20 Spark YG 593	0.6	0.8	0	1.6	-	0.7	1	0.7	0.7	1.3	1.2	0	1.5	0.5	0.2	0	1.4	0.5	0.6	0	0.6	0.3	4	2	2	3.6	5.4	2	2	6.2		6.2	4.8	6.4	5.7	4.4	3.7	2.6	3.5	5.4	3.5	1.5	0.9	0.5	0
CD337(NKp30) PE-Dazzle 594	0	5.4	0	0.4	2	0.5	0.6	0.2	2.2	2.2	3.7	0	0	1.4	4.6	0	2	1.9	0.9	0	0	0	3.9	2.5	3	6	5.4	3.3	1.9	/.4	4.9		17.9	9.7	4.1	3	1.2	0	5.3	5.7	3.8	0	0.5	0.9	0
CD4 CF594	0.5	U	0	0.6	0.8	0.6	0.5	0.5	0.3	1.5	0.7	0.7	0.9	0.3	0.5	0.7		1.6		3.2	3.6	1.5	3.3	1.9	1.8	2.8	4.3	2.1	1.7	5	6.3	5.3		6	5.9	3.3	2.3	2.1	3.7	5.5	3.5	1.7	0.6	0.5	0
CD24 PE-Alexa Fluor 610	0	1.2	1.9	1.3	1.3	1	1.4	0.9	1.8	1.4	1.9	0.3	1.4	1.6	1.7	3	3.4	2.5	1.7	3.6	0	2	5.6	3.6	4.3	7.6	9.7	3.1	2.9	7.7	10.1 1	2.6	12		7.9	4.6	3.4	2.3	8.9	10.7	7.8	3.4	1.8	2	1.8
CD95 PE-Cy5	0.2	2.5	2.2	1.2	0.7	1	1.2	0.9	1	1.7	1.3	0	1.7	0.9	0.4	1.8	2.9	2.5	1.9	0.6	1.7	0.8	3.8	4	3.6	6.5	7.9	3.4	2.9	5.2	7.3	8.2	3.5	7.3		4.6	3.2	2.7	6.3	11.5	7.4	3.7	1.8	2.3	2.2
CD25 PE-Alexa Fluor 700	0	2.4	2.3	0	0.5	0.8	1.7	1.3	0	3.1	2.8	0.4	1.9	1.1	0.2	0.8	5	3.4	2.6	1	3	1.5	2.6	1.7	3.9	6.9	11.7	5.4	3.8	3.9	4.2	7.2	2.7	5.7	3.8		4.5	4.6	5.2	9.8	5.6	4.4	2.8	2.4	2.4
CXCR3(CD183) PE-Cy7	0	2.2	2.1	0.7	0.5	0.2	0.9	1	0.9	1.1	0.9	0.5	0.5	0.4	0.1	0	1.2	2.7	3.4	0	1.4	0.5	1	0.5	1.2	2	3.5	4.7	4.5	1.7	1.7	2.2	1.1	1.8	1.3	2.1		5.8	1	2.1	1.5	1.2	1.5	2.8	2.3
HLA-DR PE-Fire 810	0.4	2.3	1.1	0	0.6	0.2	0.9	1	0	0	1.6	0.9	2.6	1.4	0.4	0	1.7	2.8	2.5	4.7	3	2.4	2.7	0.9	2	4.3	4.2	3.5	5.8	6.7	2.6	5	1.9	2.8	2.3	3	1.7		2.6	3	5.4	1.1	1.2	2.4	3.1
CD159a(NKG2A) APC	0	0	1.2	0.9	0.5	1.4	1.2	0.8	0.7	1.2	1.1	0	1.2	0.6	0.6	1.6	2.1	1.7	1.5	0	2.4	0.2	2.9	1.6	1.9	3.1	3.8	1.8	1.5	3.6	5	5.6	2.6	4.4	4.4	3.2	2.3	2		14.1	10.3	4.8	3.1	3.4	3.4
CD1c Alexa Fluor 647	1.4	0	0	0	0	0.3	1.2	0.8	1.4	4.4	1.3	1.4	2.9	1.4	0.2	1.6	0.8	1.3	0	0	10.7	1.9	2	2	0	3.3	2	0	0	5.3	7.1	7.2	3.3	5.9	5.3	2.3	2.1	3	12		21.5	6.1	3.3	5.3	4.4
CD19 Spark NIR 685	0	0	0	1.2	0	0	1.2	0.6	3.1	5.5	0	0	0	0	0	0	1.8	1.2	0.7	7.6	0	2.9	1.7	1.7	0	3.6	3.2	2.8	0.4	0	4.9	5.7	1.2	4.3	6.6	0.5	2.2	2.7	13.2	19.9		5.2	4.4	3.6	3.9
CD127 Spark Red 718	0	4.7	3.9	0	1.2	0.7	1.9	1.2	1.9	1	1.3	0.4	0	0.4	0	0	2.3	1.8	2.1	0	1.5	0	1.7	1.2	1.7	3.5	4	2	2	2.5	3.2	4.2	1.7	4.2	2.5	3.2	2.8	2.2	6.9	14.8	11.3		3.9	6.2	5.7
L/D Zombie NIR	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0		0	0
CD27 APC-H7	0	0	0.3	0.2	0	0	0.5	1.1	0	0	0	0	0	0	0	0	0	1.7	2.6	0.2	1.3	0	0	0	0	0	0.8	2	2.4	1	0.9	0.8	0	0.6	0.4	1.1	4	3.5	1.6	3.8	2.9	1.9	3.1		8.6
CD38 APC-Fire 810	0.3	0	0	0.8	0	0.4	0.5	1.3	0	0	0.5	0.1	1.2	0	0	0	0.3	1.6	2.2	0	0	0.7	0.8	0.6	0	0.9	0.3	1.8	2.4	1.7	1.8	2.1	0.6	2.1	1.7	1.2	2.4	4.4	3.4	5	4.2	2.1	1.8	6.8	

図 4.45 色のパネルの Spillover Spreading マトリックス

免疫細胞のタイプとサブタイプの分布はマーカー発現に基づいて同定され、図5に示すように階層ゲーティングを使用してそれぞれを特定した。 異なる細胞タイプに分類された分離データを図5に示します。 複雑で広がりのある多くのサブタイプの特徴付けがこのパネルで可能で あることがよくわかります。



図 5. 同定されたすべての細胞サブセットの概略図

主要な免疫細胞サブセットを同定するために使用される 手動ゲーティング戦略

PBMC は、図6に示すように、一連の染色手順に基づいたプロトコルに 従って処理されます。血小板、デブリ、ダブレット、死細胞を除外した後、 生きた CD45+ 集団内で好塩基球(1)が CD45+CD123+HLA-DR-と して分類されました。リンパ球と単球(2)は FSC-H/BSSC-H 特性に基 づいてゲーティングされました。単球(3)は CD14 と CD16 の発現に よって非古典的単球(CD14-CD16+)、中間単球(CD14+CD16+/低)、 古典的単球(CD14+CD16-)に分類されました。リンパ球ゲート(2) からは、次の集団が同定されました。CD3-TCRy δ -、CD3+TCRy δ +、 CD3+TCRyδ-(4)。CD3+TCRyδ+集団(5)はCD45RAとCCR7の 発現に基づいて特徴付けられました。CD3+TCRγδ-集団は CD3+CD56+ (NKT 様) サブセットと CD3+CD56- サブセットに類別されました(6)。 CD2 と CD8 を加えることで NKT 様細胞のさらなる分類が可能にな り、CD69 と CD57 は NKT 様細胞の早期活性化と細胞毒性を検出す るために使用されました(7)。CD3+CD56-ゲートから CD4+、CD8+、 CD4+CD8+、および CD4-CD8- T 細胞が同定されました(8)。 Treg 細 胞は CD4+ 集団から CD127 と CD25 の発現(CD127lo/-CD25hi)を 用いて同定され、これらの細胞は CD39 と CD45RA を用いてさらに分 類されました (9)。CCR7、CD45RA、CD27、CD28 により、メモリー / エフェクター CD4 および CD8 T 細胞サブセットのさらなる分類が可能に なり、CXCR3 および CCR6 により Th 集団のさらなる分類が可能になり ました。CD45RO および CD31 は CD4 T 細胞内の RTE 細胞の検出に 使用され、PD-1 および CD223 は CD4 および CD8 T 細胞の疲弊状態 の検出に使用されました(10、11)。

CD19+ および/または CD20+ 細胞 (B 細胞) は CD3-TCRvδ- 集団か ら排除されました (12)。CD19+CD20+/- 細胞はさらに lgD+CD27-、 IgD+CD27+、または IgD-CD27+/- にゲーティングされ、IgD-CD27+/ - サブセットは CD20 発現に基づいて形質芽球または lgD-メモリー B 細胞に類別され、IgD-メモリー B 細胞内で IgG および IgM 発現が 評価されました。IgD+CD27- 集団はさらに移行性 B 細胞に類別され、 IgD +CD27+ 集団はさらに MZ および IgD- メモリーのみ B 細胞に類 別されました(13)。NK 細胞は CD3-TCRγδ-HLA-DR-と定義され、初 期 NK (CD56+CD16-)、成熟NK (CD56+CD16+)、終末NK (CD56-CD16+) 細胞に分類されました(14)。樹状細胞(DC、15)は、最初 に CD3-CD19-CD56-CD14-HLA-DR+ でゲーティングすることによって 同定され、そこから CD123+ (pDC) と CD11c+DC が同定されました。 CD11c+DC はさらに CD16- と CD16+ に類別されました。次に CD1c と CD141 を使用して、CD11c+CD16- と CD11c+CD16+DC がさら に分類されました。最後に、自然リンパ球 (ILC、16) が CD3-CD19-CD20-CD14-CD123-CD127+ として同定され、CD2 および CD4 の発 現に基づいてさらに分類されました。提示されたすべてのデータは、1人 の健康なドナーの凍結 PBMC から得られたものです。



図 6. 血小板、デブリ、ダブレット、 および死細胞を除去した後に正常 なヒト PBMC サンプルで特徴付け られる細胞タイプは、単球、好塩 基球、ナチュラルキラー細胞、樹 状細胞、自然リンパ球、B リンパ 球、Tリンパ球、NKT 様リンパ球、 Treg リンパ球、および TCRγδ リ ンパ球です。分化・疲弊・活性化 の状態を評価するための他のマー カーも含まれていました。

結論

スペクトルフローサイトメトリーは、データ分解能とパネル設計の柔軟性 を向上させ、より多くのマーカーを同時に分析する機能を提供します。こ の検討で使用された Agilent NovoCyte Opteon は、使いやすい Agilent NovoExpress データ取り込みおよび解析ソフトウェアとともに、これら の利点を実現します。NovoCyte Opteon は、5 つのレーザー(UV、V、B、 Y、R)と合計 73 個の光検出器を備え、さらに高度化する実験パネル要 件を満たすことが可能で、従来のフローサイトメトリーでは見出せなかっ た分析結果を明らかにする効率的なワークフローを実現します。

参考文献

- Sen, P.; Kemppainen, E.; Oresic, M. Perspectives on Systems Modeling of Human Peripheral Blood Mononuclear Cells. *Front Mol Biosci* 2017, 4:96.
- Corkum, C.P.; Ings, D.P.; Burgess, C.; Karwowska, S.; Kroll, W.; Michalak, T.I. Immune Cell Subsets and Their Gene Expression Profiles From Human PBMC Isolated by Vacutainer Cell Preparation Tube (CPT) and Standard Density Gradient. *BMC Immunol* **2015**, 16:48.
- Kleiveland, C.R. Peripheral Blood Mononuclear Cells. In: The Impact of Food Bioactives on Health: In Vitro and Ex Vivo Models. Edited by Verhoeckx K, Coter P, Lopez-Exposito I, Kleiveland C, Lea T, Mackie A, Requena T, Swiatecka D, Wichers H. Cham (CH); 2015, 161–167.
- Lozano-Ojalvo, D.; Lopez-Fandino, R.; Lopez-Exposito, I. PBMC-Derived T Cells. In: The Impact of Food Bioactives on Health: In Vitro and Ex Vivo Models. Edited by Verhoeckx K, Coter P, Lopez-Exposito I, Kleiveland C, Lea T, Mackie A, Requena T, Swiatecka D, Wichers H. *Cham* (CH); **2015**, 169–180.
- Morath, A.; Schamel, W.W. Alphabeta and Gammadelta T Cell Receptors: Similar but Different. *J Leukoc Biol* 2020, 107(6):1045–1055.
- McBride, J.A.; Striker, R. Imbalance in the Game of T Cells: What Can the CD4/CD8 T-Cell Ratio Tell Us About HIV and Health? *PLoS Pathog* **2017**, *13*(11):e1006624.
- Golubovskaya, V.; Wu, L. Different Subsets of T Cells, Memory, Effector Functions, and CAR-T Immunotherapy. *Cancers* (Basel) **2016**, 8(3).
- 8. Sakaguchi, S.; Yamaguchi, T.; Nomura, T.; Ono, M. Regulatory T Cells and Immune Tolerance. *Cell* **2008**, *133*(5):775–787.
- 9. Raphael, I.; Nalawade, S.; Eagar, T.N.; Forsthuber, T.G. T Cell Subsets and Their Signature Cytokines in Autoimmune and Inflammatory Diseases. *Cytokine* **2015**, *74*(1):5-17.

- Pido-Lopez, J.; Imami, N.; Aspinall, R. Both Age and Gender Affect Thymic Output: More Recent Thymic Migrants in Females Than Males as They Age. *Clin Exp Immunol* **2001**, *125*(3):409–413.
- Gatinoni, L.; Speiser, D.E.; Lichterfeld, M.; Bonini, C. T Memory Stem Cells in Health and Disease. *Nat Med* **2017**, *23*(1):18– 27.
- 12. Li, H.; Tsokos, G.C. Double-Negative T Cells in Autoimmune Diseases. *Curr Opin Rheumatol* **2021**, *33*(2):163–172.
- Desfrancois, J.; Moreau-Aubry, A.; Vignard, V.; Godet, Y.; Khammari, A.; Dreno, B.; Jotereau, F.; Gervois, N. Double Positive CD4CD8 Alphabeta T Cells: A New Tumor-Reactive Population in Human Melanomas. *PLoS One* **2010**, *5*(1):e8437.
- Gonzalez-Mancera, M.S.; Bolanos, N.I.; Salamanca, M.; Orjuela, G.A.; Rodriguez, A.N.; Gonzalez, J.M. Percentages of CD4+CD8+ Double-Positive T Lymphocytes in the Peripheral Blood of Adults from a Blood Bank in Bogota, Colombia. *Turk J Haematol* **2020**, *37*(1):36–41.
- Kitchen, S.G.; Whitmire, J.K.; Jones, N.R.; Galic, Z.; Kitchen, C.M.; Ahmed, R.; Zack, J.A. The CD4 Molecule on CD8+ T Lymphocytes Directly Enhances the Immune Response to Viral and Cellular Antigens. *Proc Natl Acad Sci U S A* **2005**, *102*(10):3794–3799.
- Santegoets, S.J.; Dijkgraaf, E.M.; Bataglia, A.; Beckhove, P.; Briten, C.M.; Gallimore, A.; Godkin, A.; Goutefangeas, C.; de Gruijl, T.D.; Koenen, H.J.; et al Monitoring Regulatory T Cells in Clinical Samples: Consensus on an Essential Marker Set and Gating Strategy for Regulatory T Cell Analysis by Flow Cytometry. *Cancer Immunol Immunother* **2015**, *64*(10):1271– 1286.
- Ouaguia, L.; Morales, O.; Mrizak, D.; Ghazal, K.; Boleslawski, E.; Auriault, C.; Pancre, V.; De Launoit, Y.; Conti, F.; Delhem, N. Overexpression of Regulatory T Cells Type 1 (Tr1) Specific Markers in a Patient with HCV-Induced Hepatocellular Carcinoma. *ISRN Hepatol* **2013**, 2013:928485. <u>https://www. ncbi.nlm.nih.gov/pmc/articles/PMC4890904/</u>
- Ouaguia, L.; Mrizak, D.; Renaud, S.; Morales, O.; Delhem, N. Control of the Inflammatory Response Mechanisms Mediated by Natural and Induced Regulatory T-Cells in HCV-, HTLV-1-, and EBV-Associated Cancers. *Mediators Inflamm* 2014, 2014:564296.
- Khan, M.A.; Khan, A. Role of NKT Cells during Viral Infection and the Development of NKT Cell-Based Nanovaccines. *Vaccines* (Basel) **2021**, 9(9).

- Poli, A.; Michel, T.; Theresine, M.; Andres, E.; Hentges, F.; Zimmer, J. CD56 Bright Natural Killer (NK) Cells: an Important NK Cell Subset. *Immunology* **2009**, *126*(4):458–465.
- Melandri, D.; Zlatareva, I.; Chaleil, R.A.G.; Dart, R.J.; Chancellor, A.; Nussbaumer, O.; Polyakova, O.; Roberts, N.A.; Wesch, D.; Kabelitz, D.; et al. The γδTCR Combines Innate Immunity with Adaptive Immunity by Utilizing Spatially Distinct Regions for Agonist Selection and Antigen Responsiveness. *Nat Immunol* **2018**, *19*(12):1352–1365. <u>https://www.nature.com/articles/ s41590-018-0253-5</u>
- Yazdanifar, M.; Barbarito, G.; Bertaina, A.; Airoldi,
 I. Gammadelta T Cells: The Ideal Tool for Cancer Immunotherapy. *Cells* 2020, 9(5).
- 23. Lunt, S.Y.; Vander Heiden, M.G. Aerobic Glycolysis: Meeting the Metabolic Requirements of Cell Proliferation. *Annu Rev Cell Dev Biol* **2011**, 27:441–464.
- 24. Pieper, K.; Grimbacher, B.; Eibel, H. B-Cell Biology and Development. *J Allergy Clin Immunol* **2013**, *131*(4): 959–971.
- Jacquelot, N.; Seillet, C.; Vivier, E.; Belz, G.T. Innate Lymphoid Cells and Cancer. Nat Immunol 2022, 23(3):371–379.
- Chung, D.C.; Jacquelot, N.; Ghaedi, M.; Warner, K.; Ohashi, P.S. Innate Lymphoid Cells: Role in Immune Regulation and Cancer. *Cancers* (Basel) **2022**, *14*(9).
- Sosa Cuevas, E.; Ouaguia, L.; Mouret, S.; Charles, J.; De Fraipont, F.; Manches, O.; Valladeau-Guilemond, J.; Bendriss-Vermare, N.; Chaperot, L.; Aspord, C. BDCA1(+) cDC2s, BDCA2(+) pDCs and BDCA3(+) cDC1s Reveal Distinct Pathophysiologic Features and Impact on Clinical Outcomes in Melanoma Patients. *Clin Transl Immunology* **2020**, 9(11):e1190.
- Ouaguia, L.; Leroy, V.; Dufeu-Duchesne, T.; Durantel, D.; Decaens, T.; Hubert, M.; Valladeau-Guilemond, J.; Bendriss-Vermare, N.; Chaperot, L.; Aspord, C. Circulating and Hepatic BDCA1+, BDCA2+, and BDCA3+ Dendritic Cells Are Differentially Subverted in Patients With Chronic HBV Infection. *Front Immunol* **2019**, 10:112.

- 29. Cormican, S.; Griffin, M.D. Human Monocyte Subset Distinctions and Function: Insights From Gene Expression Analysis. *Front Immunol* **2020**, 11:1070.
- Voskamp, A.L.; Pricket, S.R.; Mackay, F.; Rolland, J.M.; O'Hehir, R.E. MHC Class II Expression in Human Basophils: Induction and Lack of Functional Significance. *PLoS One* **2013**, 8(12):e81777.
- Park, L.M.; Lannigan, J.; Jaimes, M.C. OMIP-069: Forty-Color Full Spectrum Flow Cytometry Panel for Deep Immunophenotyping of Major Cell Subsets in Human Peripheral Blood. *Cytometry A* **2020**, *97*(10):1044-1051.
- Shevchenko, Y.; Lurje, I.; Tacke, F.; Hammerich, L. Fluorochrome-Dependent Specific Changes in Spectral Profiles Using Different Compensation Beads or Primary Cells in Full Spectrum Cytometry. *Cytometry A* 2024.

www.agilent.com/lifesciences/cellanalysis

ホームページ

www.agilent.com/chem/jp

カストマコンタクトセンタ

0120-477-111 email_japan@agilent.com

本製品は一般的な実験用途での使用を想定しており、 医薬品医療機器等法に基づく登録を行っておりません。 本文書に記載の情報、説明、製品仕様等は予告なしに 変更されることがあります。

RA45406.3860069444

アジレント・テクノロジー株式会社 © Agilent Technologies, Inc. 2024 Printed in Japan, May 30, 2024 5994-7397JAJP

