



DURACLONE

RUO ANTIBODY PANELS

ADVANCED FLOW CYTOMETRY SOLUTIONS FOR CLINICAL RESEARCH

Dry, unitized, pre-formulated antibody panels for rare event detection, human adult stem cell characterization, immune function and immune system research

- Lean workflow, no antibody pipetting
- Minimized inventory management
- Dry single color reagents for compensation included*
- Ideally suited for use on the CytoFLEX and the Navios Flow Cytometers



Add whole blood sample & incubate



Lysis



Wash



Acquire



*C1101, C1102, C21857, C21858, C23406 and C75112 do not require compensation, hence do not contain compensation tubes.

DURAClone RE for Rare Event Detection

The dry DURAClone RE panels enable sensitive detection of rare abnormal cell populations in blood or bone marrow using carefully optimized and highly dosed antibody combinations.

Product - Part Number	405 nm		488 nm					633 nm		
	PB	KrO	FITC	PE	ECD	PC5.5	PC7	APC	APC-AF700	APC-AF750
RE CLB Tube - B80393 25 tests RUO	CD20	CD45	CD81	ROR-1	-	CD79b	CD19	CD5	-	CD43
RE PC Tube - B80394 25 tests RUO	CD38	CD45	CD81	CD27	-	CD19	CD200	CD138	-	CD56
RE ALB Tube - C00163 25 tests RUO	-	CD45	CD58	-	CD34	CD10	CD19	-	CD38	CD20
RE CTC Tube* - C75112 15 tests RUO	DAPI		Cyto-keratin				CD45	CD326 (EPCAM)		

*For use with PerFix CTC (C75118)

DURAClone SC for Human Adult Stem Cell Characterization

The DURAClone SC antibody panels enable rigorous human stem cell characterization in complex biological source materials as well as in enriched cellular products.

Product - Part Number	405 nm		488 nm					633 nm			
	PB	KrO	FITC	PE	ECD	PC5.5	PC7	APC	AF647	AF700	APC-AF750
SC Mesenchymal - C34369 25 tests RUO	CD31	CD14/CD19	CD90	CD73	CD34	CD146	CD105	-	-	-	CD45
SC Hematopoietic - C49589 25 tests RUO	CD90	CD45	CD38	CD49f	CD34	-	CD10	CD133	-	-	CD45RA

DURActive Stimulation Kits

The DURActive dry stimulation mixes ensure reliability of the first crucial step in cellular activation assays.

Product - Part Number	Contained Stimulating Agents	Matching DURAClone antibody panel
DURActive 1 - C11101	PMA / Ionomycin / Brefeldin A	DURAClone IF T Activation DURAClone IF T Helper Cell
DURActive 2 - C11102	PMA / Ionomycin	Detection of soluble compounds, DURAClone IF T Activation, DURAClone IF T Helper Cell (using secretion inhibitors such as Monensin)
DURActive 3 - C21857	LPS / Brefeldin A	IF Monocyte Activation

DURAClone IF for Immune Functional Cytometry

The DURAClone IF panels reveal immune function at the single cell level utilizing a streamlined and simplified staining workflow.

Product - Part Number	405 nm			488 nm				633 nm			
	PB	KrO	FITC	PE	ECD	PC5.5	PC7	APC	AF647	AF700	AF750
IF T Activation - B88649 25 tests RUO	CD4	-	IFN γ	TNF α	-	-	IL-2	-	-	CD8	CD3
IF T Helper Cell - C04666 25 tests RUO	IL-17A	-	IFN γ	-	-	-	IL-4	CD4	-	-	CD3
IF Monocytes Activation - C21858 25 tests RUO	CD14	CD45	-	HLA-DR	-	-	-	-	-	TNF α	-
IF Basophil Activation - C23406 25 tests RUO	CD63	CD45	-	CD203c	-	-	CD3	-	CD294 (CRTH2)	-	-

DURAClone IM for Immune System Research

The DURAClone IM antibody panels provide a comprehensive image of the immune system applying dry expert-proven antibody combinations that are easy to implement.

Product - Part Number	405 nm			488 nm				633 nm				
	PB	KrO	FITC	PE	ECD	PC5.5	PC7	APC	AF647	AF700	APC-AF700	APC-AF750
IM Phenotyping Basic Tube - B53309 25 tests RUO	-	CD45	CD16	CD56	CD19	-	CD14	CD4	-	CD8	-	CD3
IM B Cell Tube - B53318 25 tests RUO	IgM	CD45	IgD	CD21	CD19	-	CD27	CD24	-	-	-	CD38
IM T Cell Subsets Tube - B53328 25 tests RUO	CD57	CD45	CD45RA	CCR7	CD28	CD279 (PD-1)	CD27	CD4	-	CD8	-	CD3
IM Dendritic Cells Tube - B53351 25 tests RUO	HLA-DR	CD45	CD16	Lineage ¹	-	CD1c	CD11c	CD370 (Clec9A)	-	-	CD123	-
IM TCRs Tube - B53340 25 tests RUO	TCRV δ 2	CD45	TCR γ δ	TCR α β	HLA-DR	-	TCRV δ 1	CD4	-	CD8	-	CD3
IM Treg Tube - B53346 25 tests RUO	Helios	CD45	CD45RA	CD25	-	CD39	CD4	-	FoxP3	-	-	CD3
IM Granulocytes Tube - B88651 25 tests RUO	CD15	CD45	CD294 (CRTH2)	-	CD16	CD33	CD11b	CD274 (PD-L1)	-	-	Lineage ²	CD62L
IM Count Tube - C00162 25 tests RUO	-	-	CD45	Counting Beads	-	7-AAD	-	-	-	-	-	-
IM Innate Lymphoid Cells tube - C96081 25 tests RUO	-	CD45	CD294 (CRTH2)	Lineage ³	-	CD117	CD335 (NKp46)	CD127	-	-	-	CD161

Lineage¹ CD3 / CD19 / CD20 / CD14 / CD56

Lineage² CD3 / CD14 / CD19 / CD56

Lineage³ CD1a / CD3 / CD14 / CD16 / CD19 / CD34 / CD94 / CD123 / TCR α β / TCR γ δ / FC ϵ R1a

Keys:

PB: Pacific Blue*
KrO: Krome Orange

AF647: Alexa Fluor* 647

AF700: Alexa Fluor* 700

AF750: Alexa Fluor* 750

APC-AF700: APC-Alexa Fluor* 700

APC-AF750: APC-Alexa Fluor* 750