



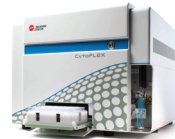
DURACLONE

RUO ANTIBODY PANELS

ADVANCED FLOW CYTOMETRY SOLUTIONS FOR CLINICAL RESEARCH

Dry, unitized, pre-formulated antibody panels for rare event detection, immune function and immune system research

- Expert-proven assays
- Lean workflow, no antibody pipetting
- Minimized inventory management
- Allowing for flexibility where needed
- 25 tests/package including 3 compensation kits
- Ideally suited for use on the CytoFlex and the Navios Flow Cytometers



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 Target cells	405 nm		488 nm				633 nm			
	PB	KrO	FITC	PE	ECD	PC5.5	PC7	APC	APC-AF700	APC-AF750
RE CLB - B80393 Abnormal CD5+ mature B cells	CD20	CD45	CD81	ROR-1	-	CD79b	CD19	CD5	-	CD43
RE PC - B80394 Abnormal plasma cells	CD38	CD45	CD81	CD27	-	CD19	CD200	CD138	-	CD56
RE ALB - C00163 Abnormal immature B cells	-	CD45	CD58	-	CD34	CD10	CD19	-	CD38	CD20

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 Pro-inflammatory T Cell Cytokines	CD4	-	IFN γ	TNF α	-	-	IL-2	-	-	CD8	CD3
IF T Helper Cell - C04666 Signature Th1/Th2/Th17 Cytokines	IL-17A	-	IFN γ	-	-	-	IL-4	CD4	-	-	CD3
IF Monocytes Activation - C21858 Innate Immune Competency	CD14	CD45	-	HLA-DR	-	-	-	-	-	TNF α	-
IF Basophil Activation - C23406 IgE-mediated Hypersensitivity	CD63	CD45	-	CD203c	-	-	CD3	-	CD294 (CRTH2)	-	-

DURAClone IM for Immune System Research

The dry 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 Leukocytes overview	-	CD45	CD16	CD56	CD19	-	CD14	CD4	-	CD8	-	CD3
IM B Cell Tube - B53318 B cell maturation	IgM	CD45	IgD	CD21	CD19	-	CD27	CD24	-	-	-	CD38
IM T Cell Subsets Tube - B53328 T cell maturation	CD57	CD45	CD45RA	CCR7	CD28	PD1	CD27	CD4	-	CD8	-	CD3
IM Dendritic Cells Tube - B53351 Subpopulations of dendritic cells	HLA-DR	CD45	CD16	Lineage ¹	-	CD11c	CD11c	Clec9A	-	-	CD123	-
IM TCRs Tube - B53340 TCR $\alpha\beta$ and $\gamma\delta$ heterodimers	TCRV δ 2	CD45	TCRV δ	TCR $\alpha\beta$	HLA-DR	-	TCRV δ 1	CD4	-	CD8	-	CD3
IM Treg Tube - B53346 Characterization of regulatory T cells	Helios	CD45	CD45RA	CD25	-	CD39	CD4	-	FoxP3	-	-	CD3
IM Granulocytes Tube - B88651 Characterization of granulocytic cells	CD15	CD45	CD294	-	CD16	CD33	CD11b	PD-L1	-	-	Lineage ²	CD62L
IM Count Tube - C00162 Viability and absolute leukocyte count	-	-	CD45	Counting Beads	-	7-AAD	-	-	-	-	-	-

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
Lineage¹: CD3 / CD19 / CD20 / CD14 / CD56
Lineage²: CD3 / CD14 / CD19 / CD56

© 2018 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries.

* Alexa Fluor and Pacific Blue are registered trademarks of Molecular Probes, Inc.