

# Cellaca PLX image cytometry system for immunophenotyping, cell counts, and viability readouts

## Multiplexity without complexity

Primary samples used by researchers in cell and gene therapy, immunotherapy, cell line development, and immuno-oncology are unique and precious due to their biological nature. The Cellaca™ PLX platform, which includes an image cytometry instrument, reagent kits, and software, is designed to increase the efficiency of sample processing, from cell staining to data readout.

The fluorescent labelled antibodies and stains in the Cellaca PLX kits are optimized for ease of use. The pre-defined assays and templates, together with the ready-to-use reagent kits, enable flow-cytometry-like data analysis with no need for titrations, extra incubations, washes, or decontamination. These features permit researchers to rapidly move processed samples to downstream applications while preserving sample integrity.



## With the Cellaca PLX platform, you can:

- Quickly perform cell purity checks and viability at the bench with Cellaca PLX immunophenotyping kits which combine CD3, CD4, and CD8 surface marker detection with viability in a single multiplex assay.
- Stain transfected cell lines with viability dyes to determine the percentage of transfected or transduced cells that are live/dead.
- Perform routine viability measurements as well as apoptosis functional assays to determine cell health.

## Cellaca PLX assay reagents and kits

1 Surface marker	Product name*	Blue	Green	Red	Far Red	Cell types	Part no. (25 Tests)	Part no. (100 Tests)
	KIRAVIA Blue 520™ anti-human CD3 Antibody*		CD3			PBMC-pure, cell lines	CS1-A0001-1	CS1-A0001-2
	Cellaca® PLX, anti-human CD3 PE Antibody			CD3		PBMC-pure, cell lines	CS1-A0002-1	CS1-A0002-2
	Cellaca® PLX, anti-human CD3 APC Antibody				CD3	PBMC-pure, cell lines	CS1-A0003-1	CS1-A0003-2
	Cellaca® PLX, anti-human CD4 PE Antibody			CD4		PBMC-pure, cell lines	CS1-A0008-1	CS1-A0008-2
	Cellaca® PLX, anti-human CD4 APC Antibody				CD4	PBMC-pure, cell lines	-	CS1-A0009-2
	Cellaca® PLX, anti-human CD8 PE Antibody			CD8		PBMC-pure, cell lines	CS1-A0012-1	CS1-A0012-2
	Cellaca® PLX, anti-human CD8 APC Antibody				CD8	PBMC-pure, cell lines	CS1-A0014-1	CS1-A0014-2
Isotype only	Product name*	Blue	Green	Red	Far Red	Cell types	Part no. (25 Tests)	Part no. (100 Tests)
	KIRAVIA Blue 520™ Mouse IgG1, κ Isotype Ctrl Antibody*		IgG1			PBMC-pure, cell lines	CS1-A0004-1	CS1-A0004-2
	Cellaca® PLX, PE Mouse IgG2a, κ Isotype Ctrl Antibody			IgG2a		PBMC-pure, cell lines	-	CS1-A0005-1
	Cellaca® PLX, APC Mouse IgG2a, κ Isotype Ctrl Antibody				IgG2a	PBMC-pure, cell lines	-	CS1-A0006-1
	Cellaca® PLX, PE Mouse IgG1, κ Isotype Ctrl Antibody			IgG1		PBMC-pure, cell lines	CS1-A0013-1**	CS1-A0013-2**
	Cellaca® PLX, APC Mouse IgG1, κ Isotype Ctrl Antibody				IgG1	PBMC-pure, cell lines	-	CS1-A0011-1
1 Surface marker + viability + isotype controls	Product name	Blue	Green	Red	Far Red	Cell types	Part no. (25 Tests)	Part no. (100 Tests)
	Cellaca® PLX, anti-human CD3 KB520 Viability Kit	Total	CD3		Dead	PBMC, PBMC-pure, cell lines	CSK-A0001-1	CSK-A0001-2
	Cellaca® PLX, anti-human CD3 PE Viability Kit	Total		CD3	Dead	PBMC, PBMC-pure, cell lines	CSK-A0002-1	CSK-A0002-2
	Cellaca® PLX, anti-human CD3 APC Viability Kit	Total	Dead		CD3	PBMC, PBMC-pure, cell lines	CSK-A0003-1	CSK-A0003-2
	Cellaca® PLX, anti-human CD4 PE Viability Kit	Total		CD4	Dead	PBMC, PBMC-pure, cell lines	CSK-A0004-1	CSK-A0004-2
	Cellaca® PLX, anti-human CD4 APC Viability Kit	Total	Dead		CD4	PBMC, PBMC-pure, cell lines	-	CSK-A0005-2
	Cellaca® PLX, anti-human CD8 PE Viability Kit	Total		CD8	Dead	PBMC, PBMC-pure, cell lines	CSK-A0006-1	CSK-A0006-2
	Cellaca® PLX, anti-human CD8 APC Viability Kit	Total	Dead		CD8	PBMC, PBMC-pure, cell lines	CSK-A0007-1	CSK-A0007-2
2 Surface markers + total + isotype controls	Product name	Blue	Green	Red	Far Red	Cell types	Part no. (25 Tests)	Part no. (100 Tests)
	Cellaca® PLX, anti-human CD3 KB520/CD4 APC Total Cell Kit	Total	CD3		CD4	PBMC, PBMC-pure, cell lines	-	CSK-A0008-2
	Cellaca® PLX, anti-human CD3 KB520/CD8 APC Total Cell Kit	Total	CD3		CD8	PBMC, PBMC-pure, cell lines	CSK-A0009-1	CSK-A0009-2
	Cellaca® PLX, anti-human CD3 APC/CD4 PE Total Cell Kit	Total		CD4	CD3	PBMC, PBMC-pure, cell lines	CSK-A0012-1	CSK-A0012-2
	Cellaca® PLX, anti-human CD3 APC/CD8 PE Total Cell Kit	Total		CD8	CD3	PBMC, PBMC-pure, cell lines	CSK-A0013-1	CSK-A0013-2
	Cellaca® PLX, anti-human CD4 PE/CD8 APC Total Cell Kit	Total		CD4	CD8	PBMC, PBMC-pure, cell lines	CSK-A0014-1	CSK-A0014-2
2 Surface markers + dead + isotype controls	Product name	Blue	Green	Red	Far Red	Cell types	Part no. (25 Tests)	Part no. (100 Tests)
	Cellaca® PLX, anti-human CD3 KB520/CD4 APC Dead Cell Kit	Dead	CD3		CD4	PBMC-pure, cell lines	-	CSK-A0016-2
	Cellaca® PLX, anti-human CD3 KB520/CD8 APC Dead Cell Kit	Dead	CD3		CD8	PBMC-pure, cell lines	CSK-A0017-1	CSK-A0017-2
	Cellaca® PLX, anti-human CD3 APC/CD4 PE Dead Cell Kit	Dead		CD4	CD3	PBMC-pure, cell lines	CSK-A0020-1	CSK-A0020-2
	Cellaca® PLX, anti-human CD3 APC/CD8 PE Dead Cell Kit	Dead		CD8	CD3	PBMC-pure, cell lines	CSK-A0021-1	CSK-A0021-2
	Cellaca® PLX, anti-human CD4 PE/CD8 APC Dead Cell Kit	Dead		CD4	CD8	PBMC-pure, cell lines	CSK-A0022-1	CSK-A0022-2
2 Surface markers + viability + isotype controls	Product name	Blue	Green	Red	Far Red	Cell types	Part no. (25 Tests)	Part no. (100 Tests)
	Cellaca® PLX, anti-human CD3 KB520/CD4 PE Viability Kit	Total	CD3	CD4	Dead	PBMC, PBMC-pure, cell lines	CSK-A0024-1	CSK-A0024-2
	Cellaca® PLX, anti-human CD3 KB520/CD8 PE Viability Kit	Total	CD3	CD8	Dead	PBMC, PBMC-pure, cell lines	CSK-A0025-1	CSK-A0025-2

## Cellaca PLX assay reagents and kits (contd).

3 Surface markers + total or dead + isotype controls	Product name	Blue	Green	Red	Far Red	Cell types	Part no. (25 Tests)	Part no. (100 Tests)
	Cellaca® PLX, anti-human CD3 KB520/CD4 PE/CD8 APC Total Cell Kit	Total	CD3	CD4	CD8	PBMC, PBMC-pure, cell lines	CSK-A0026-1	CSK-A0026-2
	Cellaca® PLX, anti-human CD3 KB520/CD4 PE/CD8 APC Dead Cell Kit	Dead	CD3	CD4	CD8	PBMC-pure, cell lines	CSK-A0027-1	CSK-A0027-2
Apoptosis detection	Product name	Blue	Green	Red	Far Red	Cell Types	Part no. (25 Tests)	Part no. (100 Tests)
	Cellaca® PLX, Caspase 3/RubyDead Apoptosis Kit	Total	Casp		Dead	Cell lines	-	CSK-A0028-2
	Cellaca® PLX, Annexin V-FITC/PI Apoptosis Kit	Total	AV	PI		Cell lines	CSK-A0029-1	CSK-A0029-2
For fluorescent proteins	Product name	Blue	Green	Red	Far Red	Cell types	Part no. (25 Tests)	Part no. (100 Tests)
	Cellaca® PLX, Hoechst/RubyDead Viability Kit	Total			Dead	Cell lines	CSK-A0030-1	CSK-A0030-2
Slides	Product name	Blue	Green	Red	Far Red	Cell types	Part no. (100 Tests)	Part no. (500 Tests)
	Cellaca® PLX, Low Fluorescence Slides					PBMC, PBMC-pure, cell lines	CHM2-ACR-001	CHM2-ACR-002

Note: The number of tests stated is based on use of the quantities described in the accompanying protocols.

\* KIRAVIA Blue™ 520 is a trademark of Sony. This product is subject to proprietary rights of Sony and is made and sold under license from Sony Corporation.

\*\*CS1-V0013-1 and CS1-V0013-2 yield a minimum of 50 and 200 tests, respectively.

