

# HTRF<sup>®</sup> SOLUTIONS FOR EACH CELL TYPE



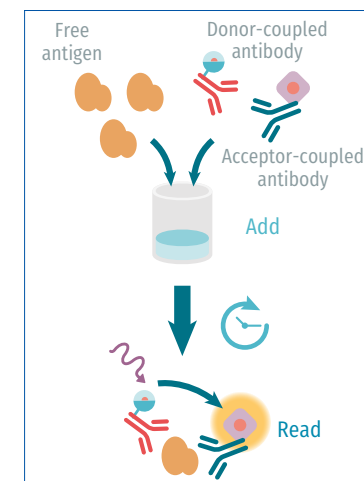
## Inflammation Cell by Cell

Inflammatory processes are complex and involve multiple cell subtypes that interact with each other. Each subtype (T cells, B cells, NK cells) expresses specific intra-cellular markers and secretes unique cytokine panels.

PerkinElmer cell-based assays offer a unique solution to monitor specific cellular activation by measuring cellular content and cytokine release.

### PerkinElmer Add-and-read Assays

- Cell-based
- Low sample volume (16  $\mu$ L)
- Intra-cellular phospho-proteins
- Intra-cellular transcription factors
- Released cytokines and biomarkers



#### ► HTRF TECHNOLOGY, A POWERFUL, EASY ALTERNATIVE TO ELISA & WESTERN BLOT

- Easy add-and-read protocol saves you time with NO washing steps
- Low sample volume consumption helps save laboratory resources

## T Cells

Subtype	Product	Size	Part
Cytotoxic CD8+	STAT1 phospho-Y701	500 TESTS	63ADK026PEG
	Granzyme B	500 TESTS	63ADK095PEG
	IFN $\gamma$	96 TESTS	62HIFNGPET
	TNF $\alpha$	96 TESTS	62HTNFAPET
T reg	STAT5 phospho-Y694	500 TESTS	64AT5PEG
	IL10	500 TESTS	62HIL10PEG
	TGF $\beta$ 1	500 TESTS	62HTGFBPEG
NKT	GM-CSF	500 TESTS	62HGMCSFPEG
	IFN $\gamma$	96 TESTS	62HIFNGPET
	TNF $\alpha$	96 TESTS	62HTNFAPET
TH1	STAT1 phospho-Y701	500 TESTS	63ADK026PEG
	IL2	96 TESTS	62HIL02PET
	IL10	500 TESTS	62HIL10PEG
	IFN $\gamma$	96 TESTS	62HIFNGPET
	TNF $\alpha$	96 TESTS	62HTNFAPET
TH2	STAT5 phospho-Y694	500 TESTS	64AT5PEG
	IL4	500 TESTS	62HIL04PEG
	IL5	500 TESTS	62HIL05PEG
	IL10	500 TESTS	62HIL10PEG
	IL31	500 TESTS	62HIL31PEG
TH9	STAT3 phospho-Y705	96 TESTS	62AT3PET
	IL10	500 TESTS	62HIL10PEG
TH17	STAT3 phospho-Y705	96 TESTS	62AT3PET
	CCL20	500 TESTS	62HCC20PEG
	IL17	96 TESTS	62HIL17PET

## B Cells

Subtype	Product	Size	Part
B-2 (Follicular)	IL2	96 TESTS	62HILPET
	IL4	500 TESTS	62HIL04PEG
	IL6	96 TESTS	62HIL06PET
	TNF $\alpha$	96 TESTS	62HTNFAPET
B reg	IL10	500 TESTS	62HIL10PEG
Be-1 (Effector 1)	IL12	500 TESTS	62HIL12PEG
	IFN $\gamma$	96 TESTS	62HIFNGPET
Be-2 (Effector 2)	IL2	96 TESTS	62HILPET
	IL4	500 TESTS	62HIL04PEG
	IL6	96 TESTS	62HIL06PET
MZ (Marginal zone)	CCL2 (MCP1)	500 TESTS	62HCCL2PEG
	CXCL10 (IP10)	500 TESTS	62HCX10PEG

## NK Cells

Product	Size	Part
AKT phospho-S473	96 TESTS	64AKSPET
ERK phospho-T202/Y204	96 TESTS	64AERPET
SLP76 phospho-S376	96 TESTS	63ADK076PET
Granzyme B	500 TESTS	63ADK095PEG
GM-CSF	500 TESTS	62HGMCSFPEG
IL10	500 TESTS	62HIL10PEG
IFN $\gamma$	96 TESTS	62HIFNGPET
TNF $\alpha$	96 TESTS	62HTNFAPET

Part numbers are for the smallest available assay format.  
Bulk sizes are available.