



# Monitor immune regulation the simple way.

## HTRF<sup>®</sup> cytokine cell-based assays

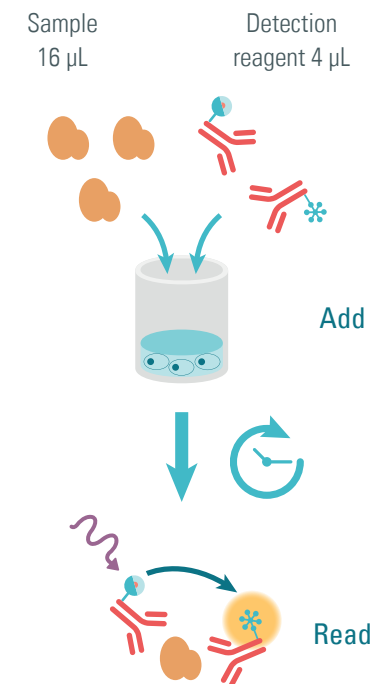
The simplest solutions always stand above the crowd.

Revvity cytokine cell-based assays offer straightforward add-and-read protocols to help you study major inflammation processes as targets or read-outs.

Now's the time to convert your ELISA testing into a truly homogeneous high throughput method. Costly, time-consuming separation and washing steps are eliminated, making this solution ideal for basic R&D, screening, and preclinical tests.

### Benefit from the most trusted TR-FRET assays to move away from ELISA:

- Small sample size
- Ideal for quantitative detection in cell lines like THP1 and primary cells such as PBMCs
- Enables multiple cytokine measurements from a single sample
- Fast and easy homogeneous format without washing steps
- Class-leading signal stability for multiple readings and reproducible results
- 96-well low volume plate included in 1x96 tests kits



## Assay specifications and ordering information

HTRF Cytokine Kits	500 Tests*	10,000 Tests*	96 Tests*
CCL2 human (MCP1)	62HCCL2PEG	62HCCL2PEH	
CCL2 mouse (MCP1)	62MCCL2PEG	62MCCL2PEI	
CCL3 human (MIP1 $\alpha$ )	62HCCL3PEG	62HCCL3PEH	
CCL4 human (MIP1 $\beta$ )	62HCCL4PEG	62HCCL4PEH	
CCL5 human (RANTES)	62HCCL5PEG	62HCCL5PEH	
CCL20 human	62HCC20PEG	62HCC20PEH	
CCL22 Human	62HCCL22PEG	62HCCL22PEH	
CXCL1 human (GRO $\alpha$ )	62HCXC1PEG	62HCXC1PEH	
CXCL10 human (IP10)	62HCX10PEG	62HCX10PEH	
G-CSF human	62HGCSFPEG	62HGCSFPEH	
GM-CSF human (CSF2)	62HGMCSFPEG	62HGMCSFPEI	
IFN alpha human	62HIFNAPEG	62HIFNAPEH	
IFN beta human	62HIFNBPEG	62HIFNBPEH	
IFN beta mouse	62MIFNBPEG	62MIFNBPEH	
IFN gamma human	62HIFNGPEG	62HIFNGPEH	62HIFNGPET**
IFN gamma mouse	62MIFNGPEG	62MIFNGPEH	
IL1 alpha human	62HIL1APEG	62HIL1APEH	
IL1 alpha mouse	62MIL1APEG	62MIL1APEH	
IL1 beta human	62HIL1BPEG	62HIL1BPEH	62HIL1BPET**
IL1 beta mouse	62MIL1BPEG	62MIL1BPEH	
IL2 human	62HIL02PEG	62HIL02PEH	62HIL02PET**
IL2 mouse	62MIL02PEG	62MIL02PEH	

HTRF Cytokine Kits	500 Tests*	10,000 Tests*	96 Tests*
IL4 human	62HIL04PEG	62HIL04PEH	
IL5 human	62HIL05PEG	62HIL05PEH	
IL6 human	62HIL06PEG	62HIL06PEH	62HIL06PET**
IL6 mouse	62MIL06PEG	62MIL06PEH	
IL8 human	62HIL08PEG	62HIL08PEH	62HIL08PET**
IL10 human	62HIL10PEG	62HIL10PEH	
IL11 human	62HIL11PEG	62HIL11PEH	
IL12 human	62HIL12PEG	62HIL12PEH	
IL12/23 p40 human	62HIL12P40PEG	62HIL12P40PEH	
IL13 human	62HIL13PEG	62HIL13PEH	
IL17 human	62HIL17PEG	62HIL17PEH	62HIL17PET**
IL17 mouse	62MIL17PEG	62MIL17PEH	
IL18 human	62HIL18PEG	62HIL18PEH	
IL22 human	62HIL22PEG	62HIL22PEH	
IL23 human	62HIL23PEG	62HIL23PEH	
IL31 human	62HIL31PEG	62HIL31PEH	
TGF beta	62HTGBPEG	62HTGBPEH	
TNF alpha cynomolgus	62CTNFAPEG	62CTNFAPEH	
TNF alpha human	62HTNFAPEG	62HTNFAPEH	62HTNFAPET**
TNF alpha mouse	62MTNFAPEG	62MTNFAPEH	

\*Based on 384-well low volume plate (20  $\mu$ L)

\*\*96-well low volume plate included

revvity