

revvity

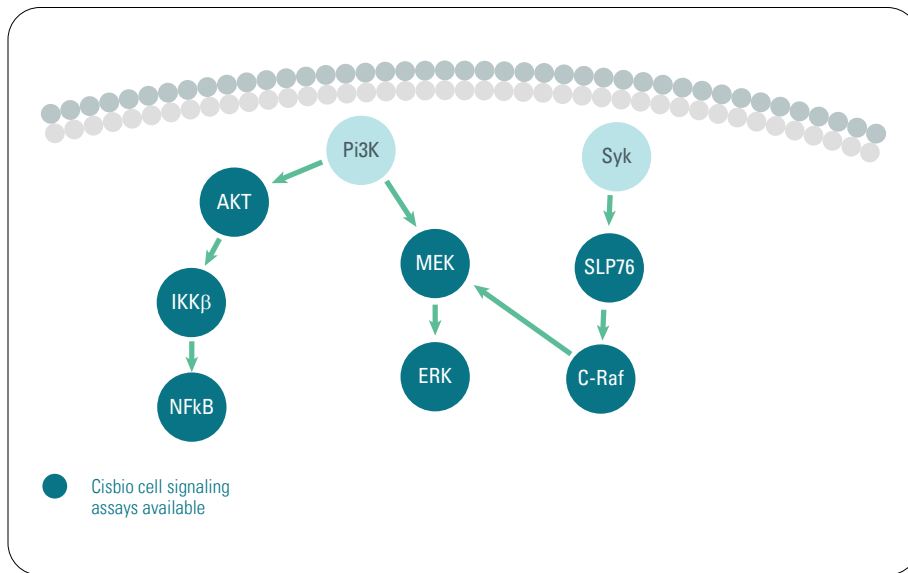
Signaling in the immune system

HTRF assays for oncology and inflammation



NK / NKT cells

Cytokines		Markers
GM-CSF (CSF2)	IL10	Granzyme B
IFN γ	TNF α	Perforin-1



Immune checkpoints

Biochemical assays to identify checkpoint blockers

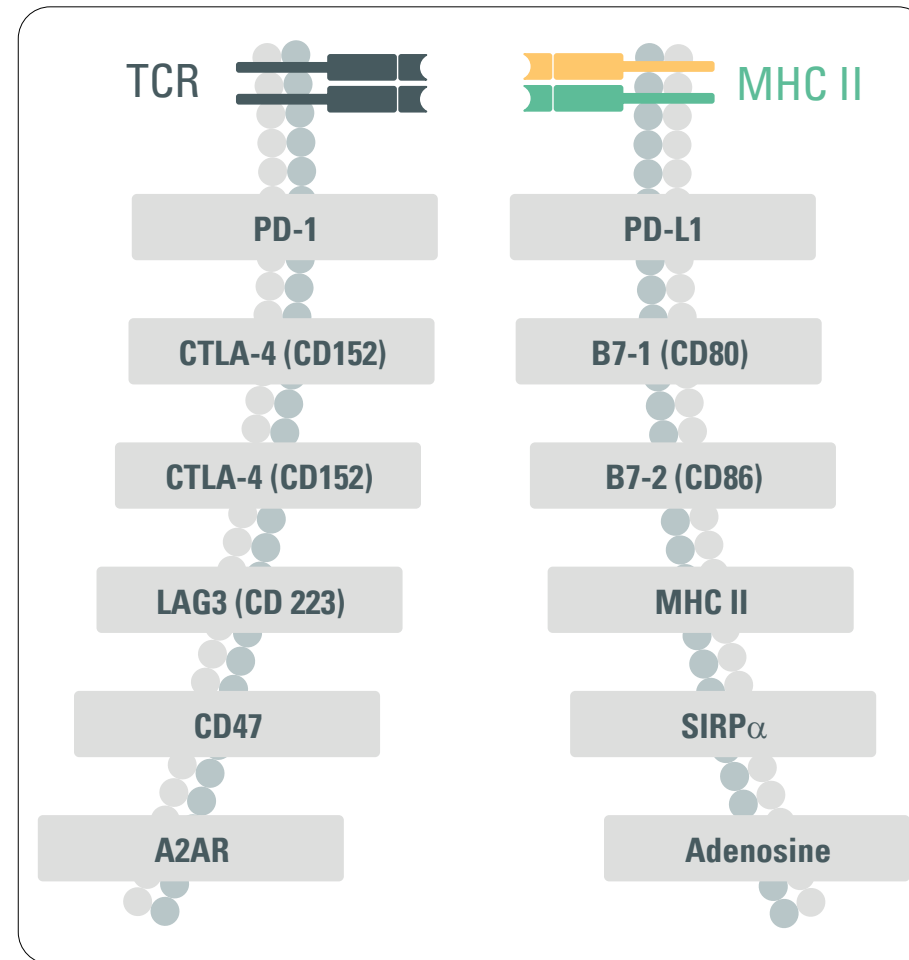


TABLE OF CONTENTS

NK / NKT cells

Immune checkpoints

T cell signaling

B cell signaling

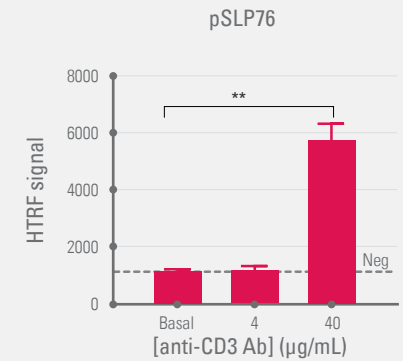
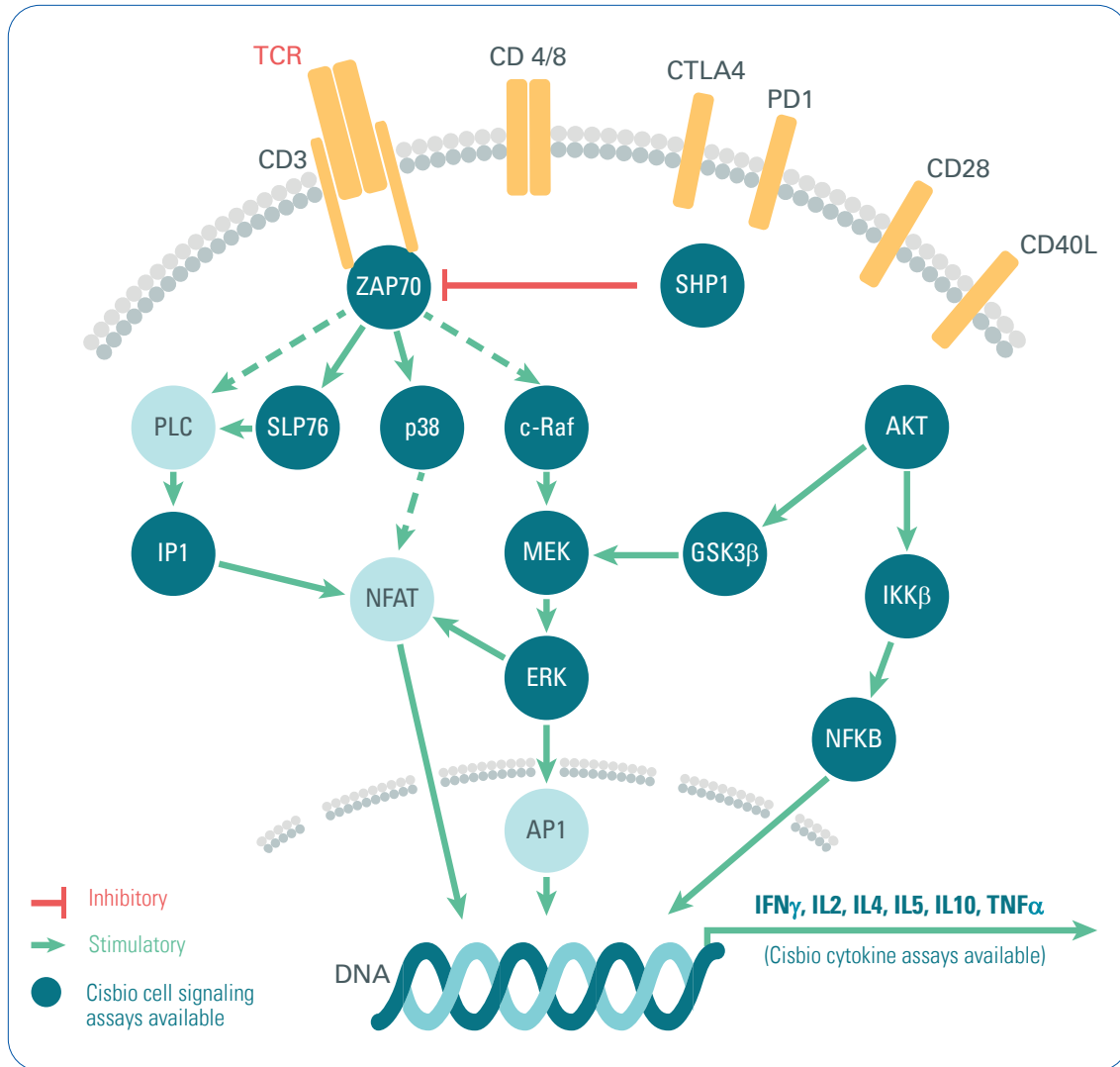
Innate immunity

Cellular phosphorylation assay

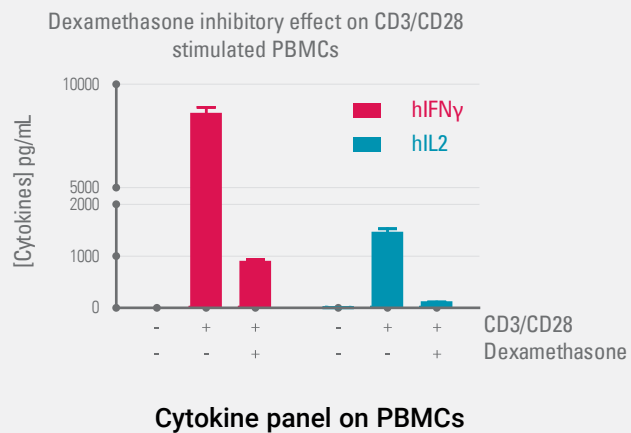
Cell-based cytokine assays

Detection

T cell signaling



Pathway monitoring in primary T cells upon CD3 stimulation
 (300,000 cells per well, 5 min CD3 stimulation)



Cytokine panel on PBMCs

TABLE OF CONTENTS

- NK / NKT cells
- Immune checkpoints
- T cell signaling
- B cell signaling
- Innate immunity
- Cellular phosphorylation assay
- Cell-based cytokine assays
- Detection



B cell signaling

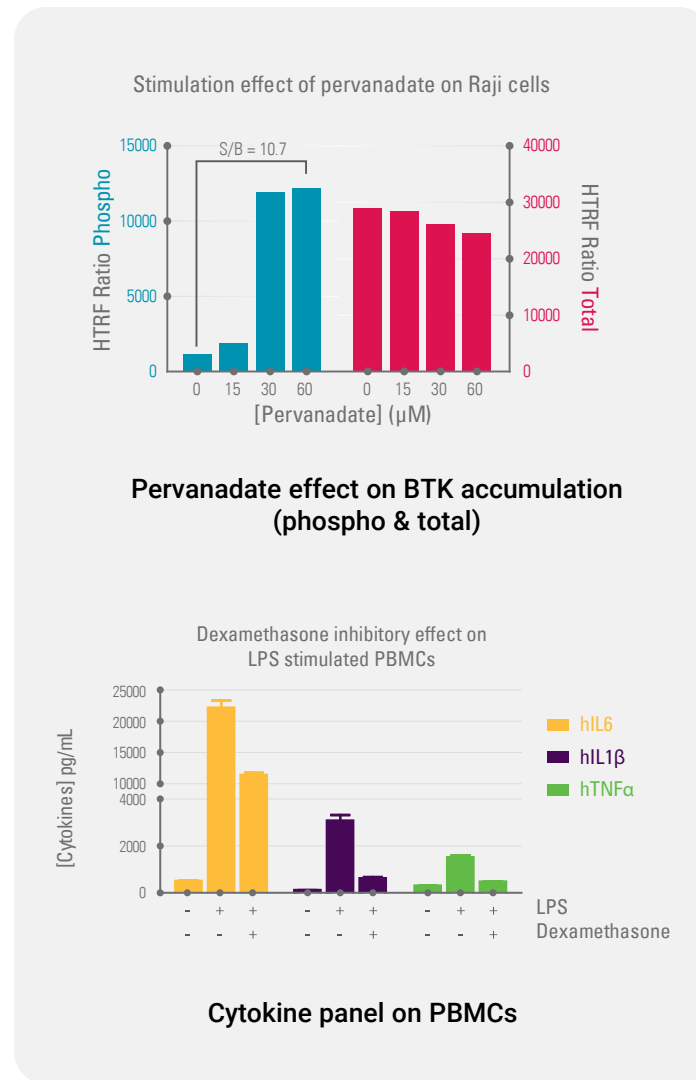
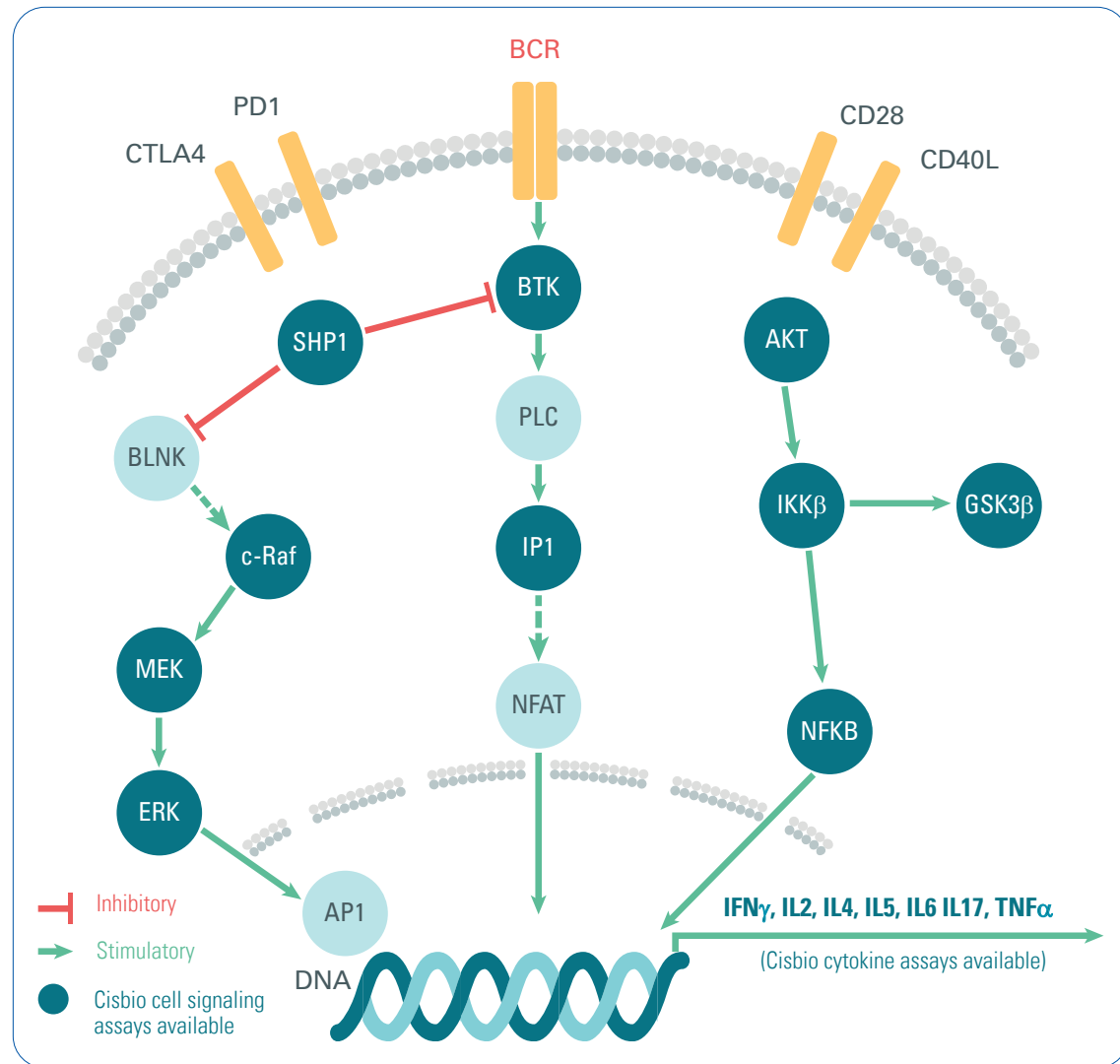


TABLE OF CONTENTS

NK / NKT cells

Immune checkpoints

T cell signaling

B cell signaling

Innate immunity

Cellular phosphorylation assay

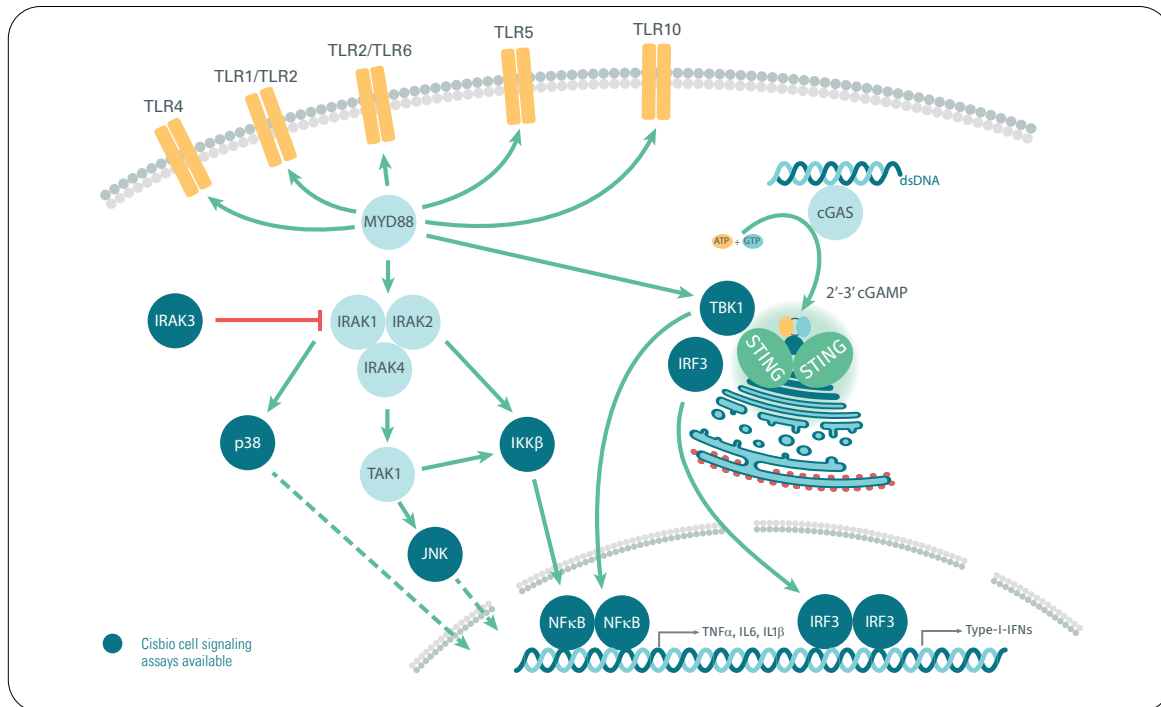
Cell-based cytokine assays

Detection



Innate immunity

Cytokines		Inflammation Markers
IL1 β	CCL3 (MIP1 α)	Histamine
IL6	CCL4 (MIP1 β)	PGE2
IL8	CCL5 (RANTES)	LTC4
IFN β	TNF α	LTB4
CXCL10 (IP-10)		



Toll-like receptor signaling pathway

STING signaling pathway

TABLE OF CONTENTS

NK / NKT cells

Immune checkpoints

T cell signaling

B cell signaling

Innate immunity

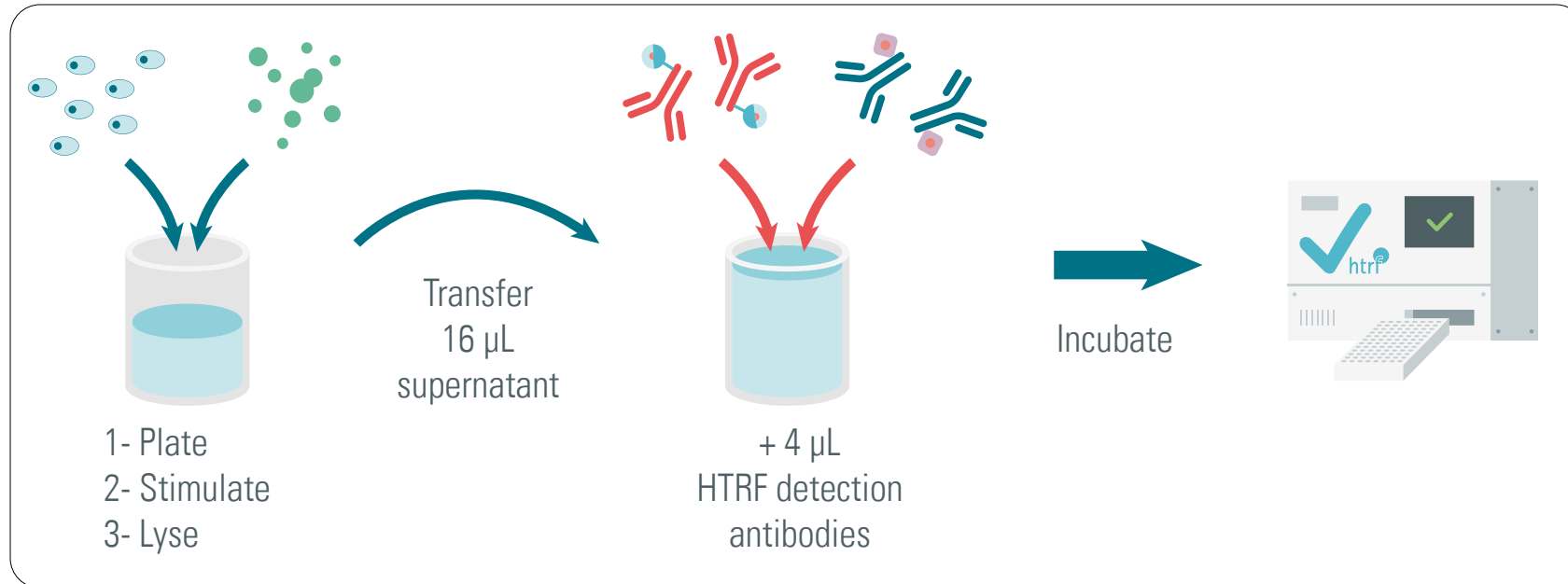
Cellular phosphorylation assay

Cell-based cytokine assays

Detection



Cellular phosphorylation assays



- Cell-based assays
- Compatible with cell lysates, tissue extracts, iPS cells, 3D cell models
- Sample volume 16 μ L
- More sensitive than Western Blot
- < 1 hr bench time
- Pharmacological studies of pathway inhibitors, agonists, and allosteric modulators

TABLE OF CONTENTS

NK / NKT cells

Immune checkpoints

T cell signaling

B cell signaling

Innate immunity

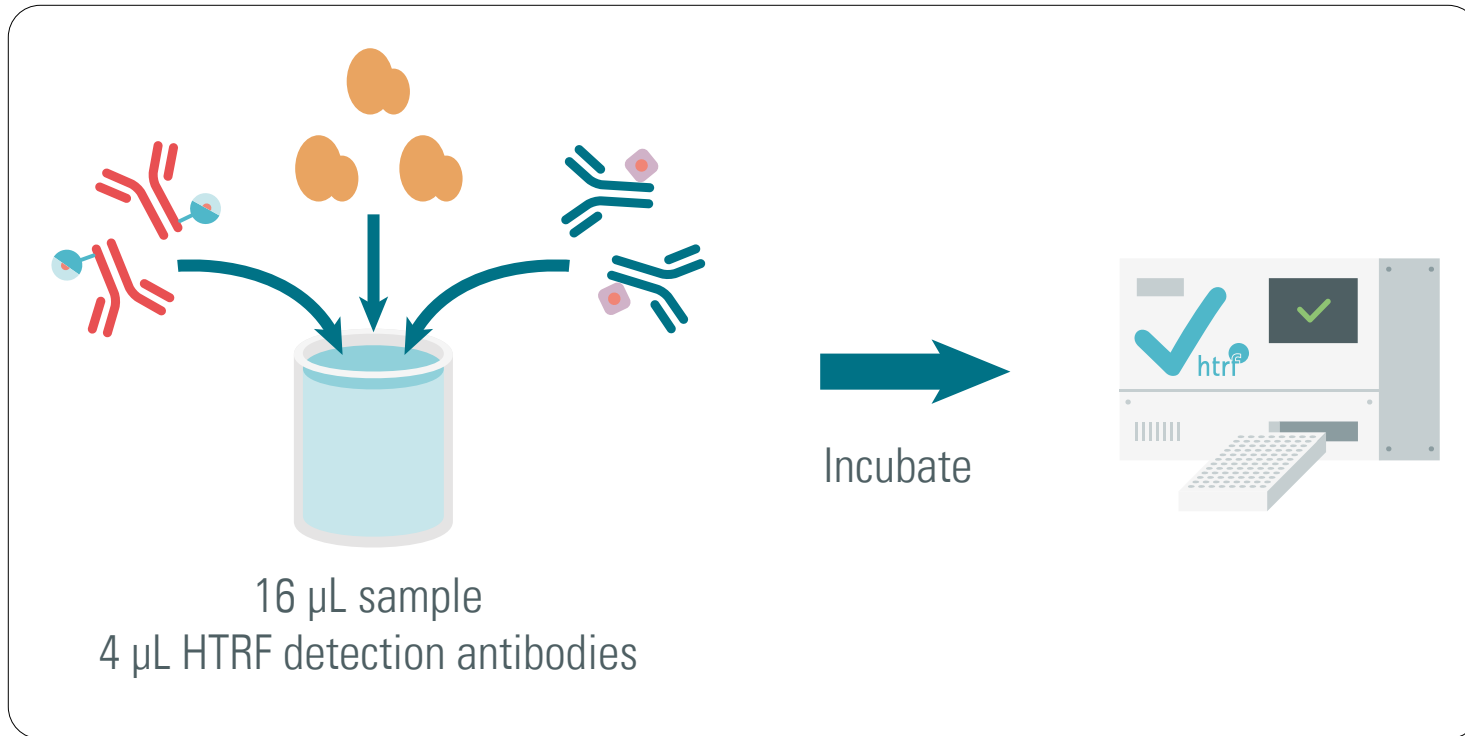
Cellular phosphorylation assay

Cell-based cytokine assays

Detection

revvity

Cell-based cytokine assays



- Sample volume 16 µL
- < 2 hrs bench time
- Validated on PBMCs and cell lines (THP1, Jurkat...)
- Compatible with Cisbio 96-well low-volume plate
- Allow multiple cytokine measurements from 1 sample

TABLE OF CONTENTS

NK / NKT cells

Immune checkpoints

T cell signaling

B cell signaling

Innate immunity

Cellular phosphorylation assay

Cell-based cytokine assays

Detection

DETECTION

Choosing the right plate reader

Choosing the suitable microplate reader ensures you'll get an optimal readout. A multimodal reader gives your research the ideal flexibility and expands the field of possibilities by measuring a wide range of technologies.

Revvity offers multimode plate readers that are certified for use with HTRF technology. Certification involves rigorous testing of plate reader with HTRF technology to ensure optimal readout. HTRF assays are also compatible with multimode readers from other providers as well.



EnVision™ Nexus™

Provides lightning speed and superior sensitivity across all established detection technologies with advanced options for ultimate performance. It is the next generation of high-throughput screening, ideal for your most demanding assays.



VICTOR™ Nivo™

Packs all the latest major detection technologies into the industry's smallest benchtop footprint. The perfect microplate reader for everyday biochemical and cell-based assays.

TABLE OF CONTENTS

NK / NKT cells

Immune checkpoints

T cell signaling

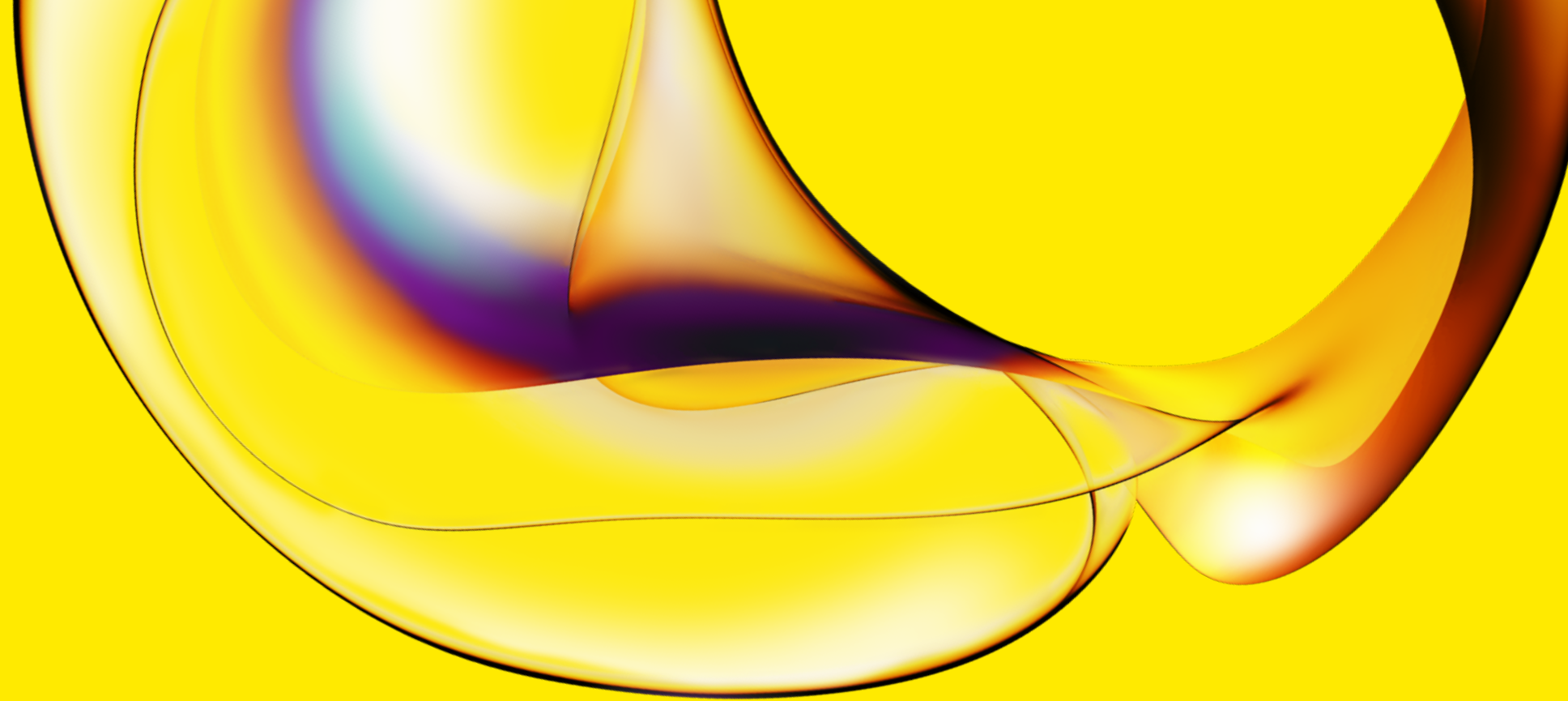
B cell signaling

Innate immunity

Cellular phosphorylation assay

Cell-based cytokine assays

Detection



www.revivity.com

revivity

Revvity, Inc.
940 Winter Street
Waltham, MA 02451 USA
www.revivity.com

For a complete listing of our global offices, visit www.revivity.com
Copyright ©2024, Revvity, Inc. All rights reserved.

1450873