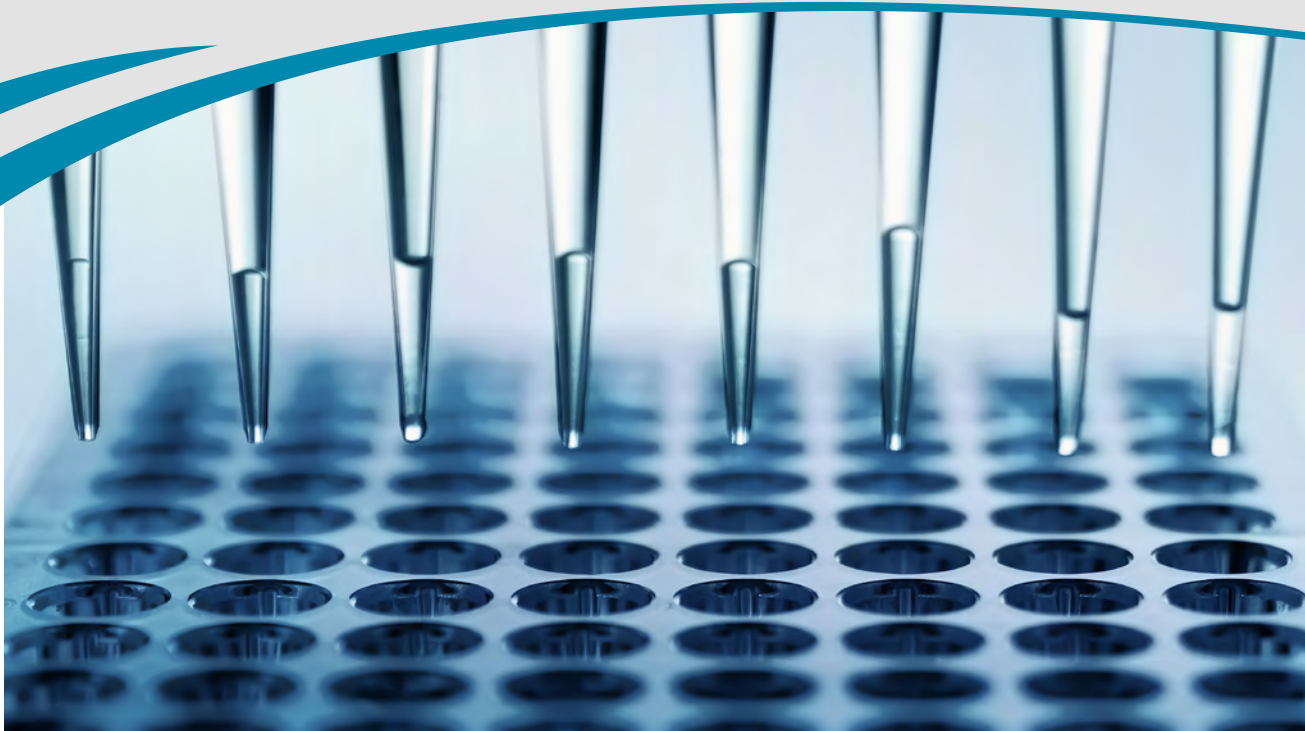


# Microplates for HTRF<sup>®</sup> immunoassays



## DESIGNED TO BE USED WITH ALL HTRF ASSAYS

Low volume and regular microplates have been carefully designed to provide the best signal detection environment possible for all HTRF assays: second messengers, phospho-proteins, cytokines, biomarkers and more....

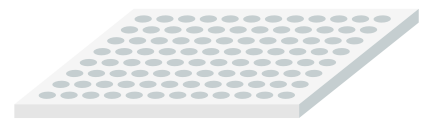
They can be used in a no-transfer protocol with suspension cells, or with a transfer step for adherent cells.

Cisbio's 96w low volume and PKI's 96w & 384w microplates can be read on all HTRF certified readers, thus making their implementation in your experiments really easy.

### At Cisbio we provide microplates that help you discover better.

- HTRF plate validated (QC with reference HTRF assays)
- Low to high volumes (10  $\mu$ L to 350  $\mu$ L)
- Can be used with all HTRF compatible readers
- For cell-based or biochemical assays

**Trying is believing!**



Color	White
Bottom	Opaque
Automation	Yes
Lid	Compatible
Sealer	Compatible (recommended)

TECHNOLOGY

## PLATE CHARACTERISTICS

	HTRF®96-well low volume plate	CulturPlate™ 96-well Microplate	CulturPlate™ 384-well Microplate	ProxiPlate™ 384 - shallow well Microplate	OptiPlate™ 384-well Microplate
Well format	96-well	96-well	384-well	384-well	384-well
Well volume	40 µL	400 µL	105 µL	28 µL	105 µL
Well shape	low volume-well	shallow-well	regular-well	regular-well	regular-well
Recommended working volume	20 µL - 30 µL	80 µL - 350 µL	24 µL - 90 µL	10 µL - 20 µL	24 µL - 90 µL
Automation Compatible	Yes	Yes	Yes	Yes	Yes
Culture	NOT tissue culture - treated	Tissue culture - treated	Tissue culture - treated	NOT tissue culture - treated	NOT tissue culture - treated
Color	White	White	White	White	White
Material	Polystyrene	Polystyrene	Polystyrene	Polystyrene	Polystyrene
Sterile	No	Yes	Yes	No	No
Provider	Cisbio	PerkinElmer	PerkinElmer	PerkinElmer	PerkinElmer

PART NUMBER#	PRODUCT DESCRIPTION	NUMBER OF PLATES#
66PL96001	HTRF 96 well low volume plate	1 plate
66PL96005	HTRF 96 well low volume plate	5 plates
66PL96025	HTRF 96 well low volume plate	25 plates
6005680	CulturPlate-96, White Opaque 96-well Microplate	50 plates
6005688	CulturPlate-96, White Opaque 96-well Microplate	160 plates
6007680	CulturPlate-384, White Opaque 384-well Microplate	50 plates
6007688	CulturPlate-384, White Opaque 384-well Microplate	160 plates
6008280	ProxiPlate-384 Plus, White 384-shallow well Microplate	50 plates
6008289	ProxiPlate-384 Plus, White 384-shallow well Microplate	200 plates
6007290	OptiPlate-384, White Opaque 384-well Microplate	50 plates
6007299	OptiPlate-384, White Opaque 384-well Microplate	50 plates

PART NUMBER#	PRODUCT DESCRIPTION	NUMBER OF TOPSEALS
6050185	TopSeal-A PLUS	box 100 items

### FOR MORE INFORMATION

Europe and other countries +33(0)466 796 705 U.S. and Canada 1-888-963-4567

China +86 10 8080 9288 Japan +81-(0)43-306-8712

Visit [www.cisbio.com](http://www.cisbio.com) to find a list of our regional distributors



cisbio