

DESIGNED TO BE
USED WITH
ALL HTRF ASSAYS



Microplates for HTRF® Immunoassays

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.

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



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

At PerkinElmer we Provide Microplates That Help You Discover Better.

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

Trying is Believing!

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	PerkinElmer	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	200 plates

Part Number#	Product Description	Number of Topseals
6050185	TopSeal-A PLUS	Box 100 items

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
P: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright ©2022, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.