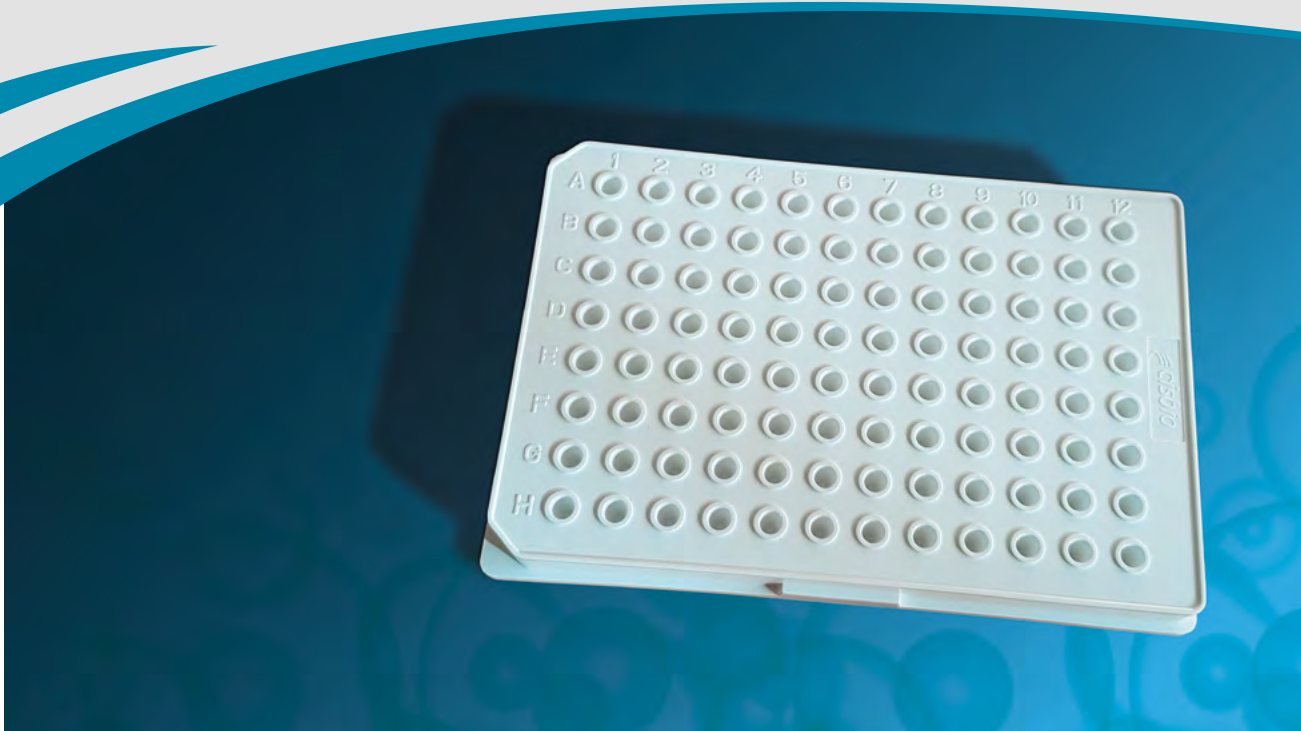


# HTRF® 96-well low volume plate



## ALL THE BENEFITS OF HTRF WITH THE EASE OF USE OF AN ELISA PLATE!

We have designed these plates to simplify your access to HTRF by enabling easy pipetting and handling while keeping the advantages of the HTRF technology:

- HTRF validated (QC with reference HTRF assays)
- Low volume (20µl)
- Compatible with all HTRF readers
- For cell-based or biochemical assays

**Trying is believing**

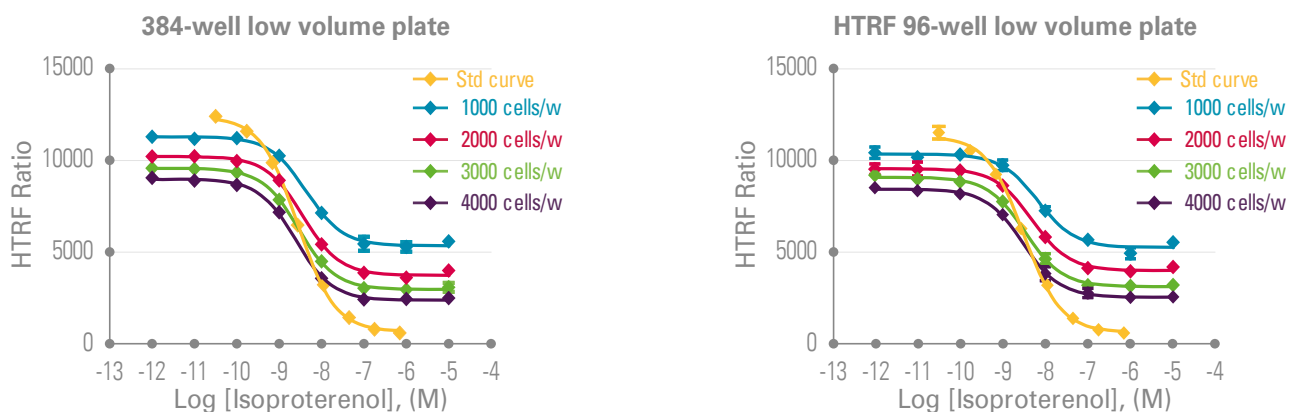


## PLATE CHARACTERISTICS

Format	96 well
Color	White
Bottom	Opaque
Maximum volume / well	40 $\mu$ L
Recommended working volume / well	20-30 $\mu$ L
Culture	NOT culture treated
Lid	Compatible
Sealer	Compatible (and recommended)
Reader settings	96 half-area (except Envision)
Validation	Inter-batch & intra-batch

## PLATE VALIDATION

5  $\mu$ L of HEK293 cells expressing the  $\beta$ 2 adrenergic receptor were dispensed side-by-side in the HTRF 96-well low volume plate and in a 384-well low volume plate. Cells were stimulated for 45 min at 37°C with 5  $\mu$ L of isoproterenol agonist, then 5  $\mu$ L of cAMP-d2 and 5  $\mu$ L of anti cAMP- cryptate from the HTRF cAMP Gs Dynamic kit were added. The HTRF signal was then recorded after a 1 hour incubation at RT. (Fig.1)



	S/B			
Cells/w	1000	2000	3000	4000
384-well low volume plate	2.1	2.8	3.2	3.6
HTRF 96-well low volume plate	2.1	2.5	2.9	3.4

	EC <sub>50</sub> (nM)			
Cells/w	1000	2000	3000	4000
384-well low volume plate	4.4	3.6	2.9	2.4
HTRF 96-well low volume plate	6.7	4.8	3.3	3.0

Fig.1 - Compound potency, EC50 (nM)

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
66PL96100	HTRF 96 well low volume plate	100 PLATES

## INTERACTION IS EVERYTHING



From the chemistry of our assays to the collaborative nature of our customer care relationships, Cisbio's approach is based on the belief that "interaction is everything". Talk to us and discover a partner dedicated to providing exactly what you need:

- Convenient ready-to-use kits
- Custom assay capabilities
- Trusted and proven HTRF technology
- Complete confidentiality
- Expert technical support

## FOR MORE INFORMATION

Europe and other countries +33(0)466 796 705 U.S. and Canada 1-888-963-4567 China +86 21 5018 9880 Japan +81-(0)43-306-8712 Visit [Cisbio.com](http://Cisbio.com) to find a list of our regional distributors

