

## HTRF® Research Reagents

## HTRF Bind. Domain kit C - 10K pts

Product Number : 62BDCPEH Lot Number : 06C

Expiration date : 2026-01-06 Document version : 1

Reagents :	Lot n°
HTRF Bind. Domain - Diluent Buf.- 200 mL	12RA
HTRF Bind. Domain - Detect. Buf.2-130 mL	14RA
HTRF Bind. Domain kit - 10K pts Tb Cryptate antibody	03RA
HTRF Bind. Domain kit - 10K pts XL665 reagent	06RA

## Results :

Incubation :

Plate :

Reader : Pherastar Plus

Fitting model :

## Assay characteristics :

[pept 001-  
biot]en nMf

	strepta-XL665 ref lot 06A		
	GST-K Tb réf lot 03A		
	Moy Ratio	CV en %	Moy Ratio Spé
Blanc	822	2%	-5
0.03	830	1%	18
0.10	843	1%	3
0.30	940	2%	119
0.91	1080	0%	245
2.74	2374	7%	1545
8.23	5877	3%	5040
24.69	11046	#DIV/0!	10166
74.07	28021	7%	27095
222.22	37352	0%	36291
666.67	46559	3%	45196
2000.00	65343	5%	62808

HTRF Bind. Domain kit C - 10K pts - Lot n° : 06C

☒ Accepted☐ Rejected

Quality control date : 2024-06-10 Supervisor's initials : ML

The information provided in this document is valid for the specified lot number and date of analysis. This information is for reference purposes only and does not constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume any liability for any errors or damages arising from the use of this document or the product described herein. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.