

HTRF® Research Reagents

HTRF Bind. Domain kit A - 500 pts

Product Number : 62BDAPEG Lot Number : 02F

Expiration date : 27/09/2024 Document version : 1

Reagents :		Lot n°
HTRF Bind. Domain - Diluent buf.- 20 mL		06RD
HTRF Bind. Domain - Detect. Buf.1- 20 mL		11RA
HTRF Bind. Domain kit - 500 pts d2 reagent		02RA
HTRF Bind. Domain kit - 500 pts Eu Cryptate antibody		02RA

Results :

Incubation :

Plate :

Reader : Pherastar Plus

Fitting model :

Assay characteristics :

[pept 001-biot]en nMf	test strepta-d2 lot 02		
	GST-K lot 02		
	Moy ratio	CV en %	moy ratio Spé
Blanc	344	2%	-18
0,03	371	2%	21
0,10	394	1%	36
0,30	446	6%	69
0,91	691	#DIV/0!	331
2,74	1207	4%	840
8,23	2715	5%	2337
24,69	6304	1%	5901
74,07	11584	1%	11102
222,22	16894	1%	16166
666,67	20468	2%	19108
2000,00	24966	1%	21666

HTRF Bind. Domain kit A - 500 pts - Lot n° : 02F

☒ Accepted☐ Rejected

Quality control date : 2022-10-06 Supervisor's initials : AC

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.