



HTRF® Research Reagents

HTRF Bind. Domain kit B - 10K pts

Product Number :62BDBPEH

Lot Number :10B

Expiration date :2025-02-15

Document version :1

Reagents :	Lot n°
HTRF Bind. Domain - Diluent Buf.- 200 mL	12RA
HTRF Bind. Domain - Detect. Buf.1-130 mL	13RA
HTRF Bind. Domain kit - 10K pts Eu Cryptate antibody	07RA
HTRF Bind. Domain kit - 10K pts XL665 reagent	10RA

Results :
Incubation : Overnight at RT
Plate : 384 wells small volume white plate 20 µL
Reader : /
Fitting model : /

Assay characteristics :

gamme nég		BRD2 (1)		
strepta-XL ref lot 10A				
GST-K lot 07A				
Moy ratio	CV en %	Moy ratio	CV en %	moy ratio Spé
369	2%	351	5%	-18
362	3%	378	2%	15
357	4%	395	4%	38
357	2%	470	7%	113
370	2%	707	8%	337
367	2%	1577	2%	1210
386	6%	5503	#DIV/0!	5117
404	6%	10504	2%	10100
449	1%	21929	0%	21480
608	3%	27860	1%	27252
1022	5%	34871	2%	33849
2520	1%	52207	1%	49687

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

☒ Accepted

☐ Rejected

Quality control date :2024-04-29

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.

