

## HTRF® Research Reagents

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

Product Number : 62BDBPEH Lot Number : 10C

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

| Reagents :                                           |  | Lot n° |
|------------------------------------------------------|--|--------|
| HTRF Bind. Domain - Diluent Buf.- 200 mL             |  | 12RA   |
| HTRF Bind. Domain - Detect. Buf.1-130 mL             |  | 13B    |
| 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      |         | 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 : 2025-03-07 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.