

HTRF® Research Reagents

HTRF (h/m) RBM39 Total Kit - 500 pts

Product Number : 64RBM39TPEG Lot Number : 01D

Expiration date : 2026-04-03 Document version : 1

| Reagents : | Lot n° |
|---|--------|
| HTRF (h/m) RBM39 Total Kit - 500 pts d2 antibody | 01B |
| HTRF (h/m) RBM39 Total Kit - 500 pts Eu Cryptate antibody | 01B |
| HTRF (h/m) RBM39 Total Kit - Ctrl Lysate | 02A |
| HTRF P-T prot. - Lysis Buf.4 (4X) 2 mL | 14B |
| HTRF P-T prot. - Block. reag.(100X) 0.3 mL | 28B |
| HTRF P-T prot. - Detect. Buf. 2 mL | 32B |

Results :

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

Assay characteristics :

| | Found value | Specifications Rev 01 |
|----------------|-------------|-----------------------|
| Negative | 491 | / |
| Control Lysate | 3975 | / |
| S/B | 8.1 | >5 |
| | / | / |

HTRF (h/m) RBM39 Total Kit - 500 pts - Lot n° : 01D

 Accepted

 Rejected

| | |
|-----------------------------------|----------------------------|
| Quality control date : 2024-08-01 | 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.