

HTRF® Research Reagents

HTRF Hu. Amyloid b1-40 kit - 500 pts

Product Number : 62B40PEG Lot Number : 11F

Expiration date : 2025-06-06 Document version : 1

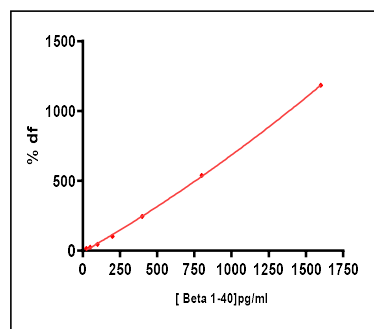
| Reagents : | Lot n° |
|---|--------|
| HTRF Diluent 1 - 20 mL | 20RA |
| HTRF Detect. Buf. 2 - 7 mL | 16RA |
| HTRF Hu. Amyloid b1-40 kit - Std | 09RA |
| HTRF Hu. Amyloid b1-40 kit - 500 pts Eu Cryptate antibody | 11RA |
| HTRF Hu. Amyloid b1-40 kit - 500 pts XL665 antibody | 11RA |

Results :

Incubation : Overnight at 4°C
 Plate : 384 wells small volume white plate 20 µL
 Reader : /
 Fitting model : /

Assay characteristics :

| B1-40 | Conc. (pg/ml) | DF % | DF Accepted |
|-------|---------------|-------|-------------|
| Std 1 | 25 | 16% | 10-25 |
| Std 2 | 50 | 24% | 20-43 |
| Std 3 | 100 | 46% | 40-74 |
| Std 4 | 200 | 102% | 83-146 |
| Std 5 | 400 | 245% | 171-299 |
| Std 6 | 800 | 539% | 374-657 |
| Std 7 | 1600 | 1184% | 837-1388 |



| | Found value | Specifications |
|------------------|-------------|----------------|
| Detection limite | 6,2 | <=10 pg/ml |

HTRF Hu. Amyloid b1-40 kit - 500 pts - Lot n° : 11F

Accepted

Rejected

| | |
|-----------------------------------|----------------------------|
| Quality control date : 2023-04-24 | 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.