

HTRF® Research Reagents

HTRF Hu. Amyloid b1-42 kit - 500 pts

Product Number : 62B42PEG Lot Number : 04D

Expiration date : 2025-03-23 Document version : 1

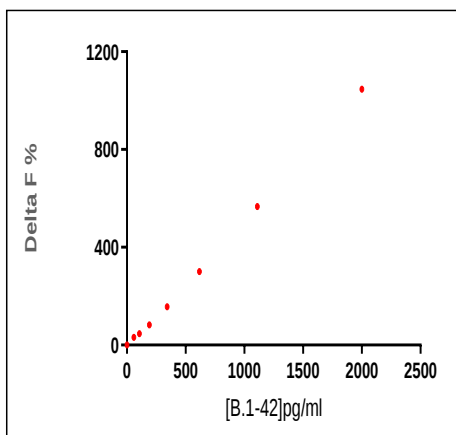
Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	21RA
HTRF Detect. Buf. 3 - 1.5 mL	30AR
HTRF Hu. Amyloid b1-42 kit - Std	01RA
HTRF Hu. Amyloid b1-42 kit - 500 pts d2 antibody	04RA
HTRF Hu. Amyloid b1-42 kit - 500 pts Eu Cryptate antibody	04RA

Results :

Incubation : Overnight at RT
Plate : 384 wells small volume white plate 20 µL
Reader : Pherastar Plus
Fitting model : 4 Parameter Logistic (4PL)

Assay characteristics :

	Conc (pg/ml)	Delta F %	Delta F % Accepted Range
Std 0	0	0%	0-0
Std 1	59	37	25-37
Std 2	106	67	45-67
Std 3	191	104	78-118
Std 4	343	169	136-204
Std 5	617	288	245-368
Std 6	1111	472	434-651
Std 7	2000	775	754-1132



HTRF Hu. Amyloid b1-42 kit - 500 pts - Lot n° : 04D

☒ Accepted☐ Rejected

Quality control date : 2023-11-30 Supervisor's initials : FMV

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.