

HTRF® Research Reagents

HTRF Hu. Amyloid b1-40 kit - 500 pts

Product Number : 62B40PEG Lot Number : 11G

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

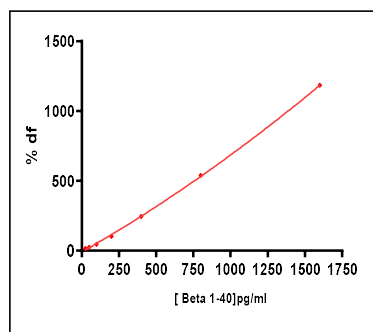
Reagents :	Lot n°
HTRF Diluent 1 - 20 mL	20RA
HTRF Detect. Buf. 2 - 7 mL	17AR
HTRF Hu. Amyloid b1-40 kit - Std	09RA
HTRF Hu. Amyloid b1-40 kit - 500 pts Eu Cryptate antibody	11RA
HTRF Hu. Amyloid b1-40 kit - 500 pts XL665 antibody	11RA

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° : 11G

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.