

HTRF® Research Reagents

HTRF BDNF kit - 500 pts

Product Number : 62BDNFPEG Lot Number : 02B

Expiration date : 2024-09-01 Document version : 1

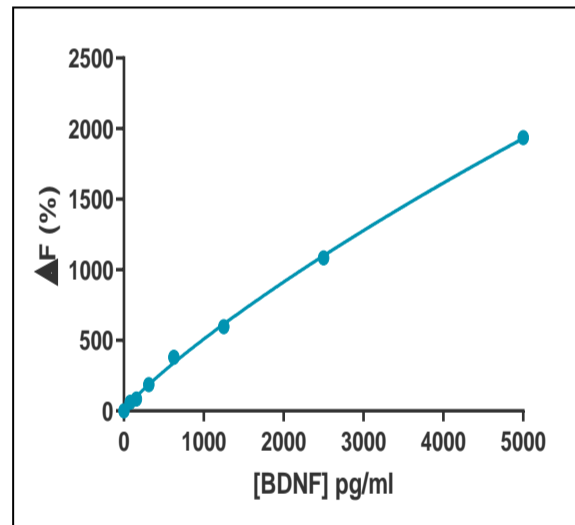
Reagents :	Lot n°
HTRF Diluent 5 (5X) - 2 mL	21RA
HTRF Detect. Buf. 3 - 1.5 mL	30AR
HTRF BDNF kit - Std	01RA
HTRF BDNF kit - 500 pts Eu Cryptate antibody	01RA
HTRF BDNF kit - 500 pts XL665 antibody	02AR

Results :

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

Assay characteristics :

BDNF	Conc. (pg/ml)	DF %	DF Accepted Range Provisoire
Std 1	78,0	37%	34-51
Std 2	156,0	88%	70-105
Std 3	313,0	142%	141-211
Std 4	625,0	289%	275-413
Std 5	1250,0	577%	513-769
Std 6	2500,0	1127%	942-1413
Std 7	5000,0	1919%	1529-2294



	Found value	Specifications Rev 01
Detection limite	10	≤ 78 pg/ml

HTRF BDNF kit - 500 pts - Lot n° : 02B

Accepted

Rejected

Quality control date : 2024-01-29	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.