

HTRF® Research Reagents

HTRF (h/m) Ataxin 2 Kit - 500 pts

Product Number : 64ATA2PEG Lot Number : 02B

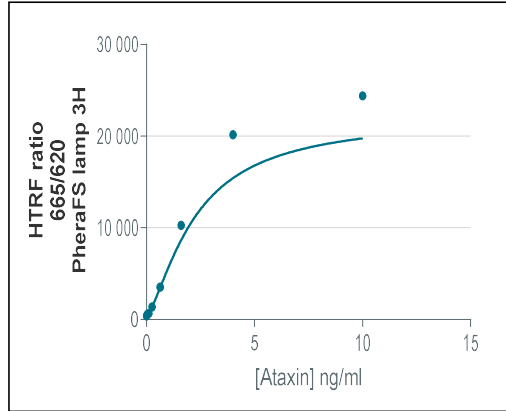
Expiration date : 2026-03-20 Document version : 1

Reagents :	Lot n°
HTRF Detect. Buf. 3 - 1.5 mL	30B
HTRF (h/m) Ataxin 2 Kit - Std	01A
HTRF (h/m) Ataxin 2 Kit - 500 pts d2 antibody	02A
HTRF (h/m) Ataxin 2 Kit - 500 pts Eu Cryptate antibody	01B
HTRF P-T prot. - Lysis Buf.1 (4X) 2 mL	35B
/	/

Results :

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

	Conc (ng/ml)	Ratio	Ratio Accepted Range
Std 0	0	391	291-482
Std 1	0.041	501	341-658
Std 2	0.1	648	488-864
Std 3	0.256	1350	928-1897
Std 4	0.64	3516	2448-4138
Std 5	1.6	10266	5807-12446
Std 6	4	20146	11340-22599
Std 7	10	24383	12868-25444



HTRF (h/m) Ataxin 2 Kit - 500 pts - Lot n° : 02B

Accepted

Rejected

Quality control date : 2024-02-15	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.