

HTRF® Research Reagents

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

Product Number : 64ATA2PEG Lot Number : 01A

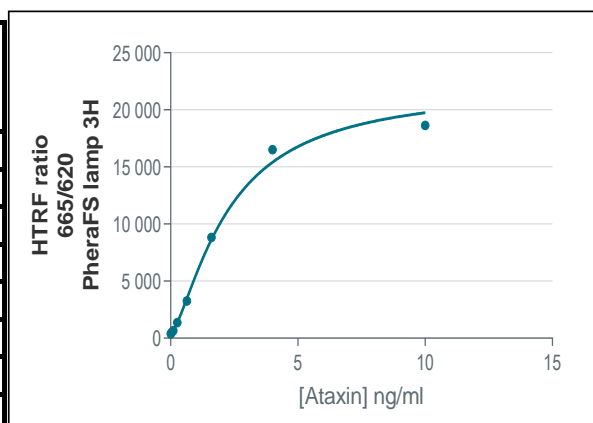
Expiration date : 12/07/2023 Document version : 1

Reagents :	Lot n°
HTRF (h/m) Ataxin 2 kit - 500 pts d2 antibody	01A
HTRF (h/m) Ataxin 2 kit - 500 pts Eu cryptate antibody	01A
HTRF (h/m) Ataxin 2 kit - Std	01A
HTRF P-T prot. Lysis Buffer 1 (4X) 2 mL	33A
HTRF Detect Buf 3 -1,5mL	29A
HTRF (h/m) Ataxin 2 kit - modop	1

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	383	291-482
Std 1	0,041	487	341-658
Std 2	0,1	665	488-864
Std 3	0,256	1370	928-1897
Std 4	0,64	3254	2448-4138
Std 5	1,6	8815	5807-12446
Std 6	4	16503	11340-22599
Std 7	10	18644	12868-25444



HTRF (h/m) Ataxin 2 kit - 500 pts - Lot n° : 01A

 Accepted

 Rejected

Quality control date : 2023-09-29

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.