

HTRF® Research Reagents

HTRF FcRn Bind. kit - 100 pts

Product Number : 64FCRNPET Lot Number : 08D

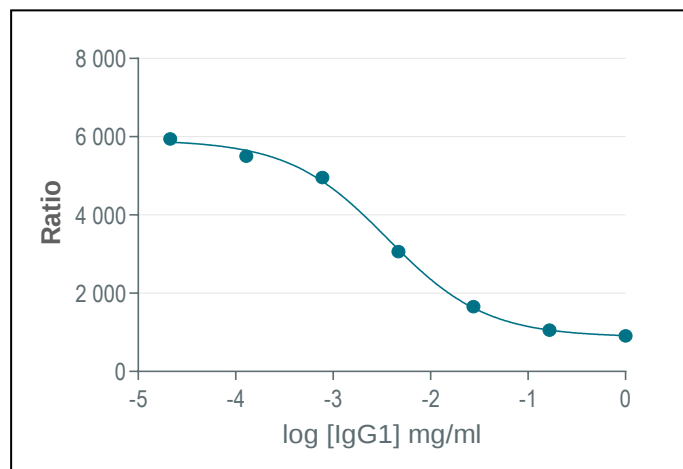
Expiration date : 2026-02-27 Document version : 1

Reagents :	Lot n°
HTRF Diluent 10 - 10 mL	05AR
HTRF Detect. Buf. 13 - 1.2 mL	05AR
HTRF FcRn Bind. kit - Std	06AR
HTRF FcRn Bind. kit - 100 pts d2 antibody	08A
HTRF FcRn Bind. kit - 100 pts Tb Cryptate reagent	08A

Results :

Incubation : 1h at RT
Plate : 384 wells small volume white plate 20 µL
Reader : Pherastar Plus
Fitting model : 4 Parameter Logistic (4PL)

	Conc (µg/ml)	ratio	ratio Accepted Range
Std 0	0	6079	5962-8554
Std 1	0,02	5 946	5208-7945
Std 2	0,13	5 502	4394-7603
Std 3	0,77	4 954	3159-6854
Std 4	5	3 066	2379-4266
Std 5	28	1 655	1420-1926
Std 6	167	1 058	1003-1171
Std 7	1000	908	738-1075



Assay characteristics :

	Found value	Specifications Rev 01
Ratio max	6079	5962-8554
EC50	3,6	1-5 µg/mL

HTRF FcRn Bind. kit - 100 pts - Lot n° : 08D

☒ Accepted

☐ Rejected

Quality control date : 2025-01-20 Supervisor's initials : PL

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.