

HTRF® Research Reagents

HTRF Estradiol kit - 500 pts

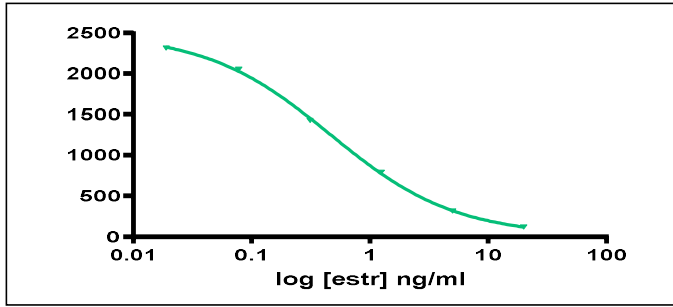
Product Number : 62ESTPEG Lot Number : 04Q

Expiration date : 2026-01-11 Document version : 1

Reagents :	Lot n°
HTRF Diluent 2 - 20 mL	10RA
HTRF Detect. Buf. 2 - 7 mL	17AR
HTRF Estradiol kit - Std	08AR
HTRF Estradiol kit - 500 pts Eu Cryptate antibody	04RD
HTRF Estradiol kit - 500 pts XL665 reagent	04RD

Results :
 Incubation : 2h at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : Pherastar Plus
 Fitting model : nonlinear regression, sigmoidal dose response (variable slope)

Estradiol	Conc. (ng/ml) _i	DF
Std 0	0	2670%
Std 1	0.019	2356%
Std 2	0.078	2019%
Std 3	0.312	1485%
Std 4	1.25	780%
Std 5	5	308%
Std 6	20	107%



Assay characteristics :

	Found value	Specifications Rev 01
Detection limit	0.006	< 0.019 ng/ml
Negative	2670	>1250%

HTRF Estradiol kit - 500 pts - Lot n° : 04Q

Accepted

Rejected

Quality control date :	2024-03-08	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.