

HTRF® Research Reagents

HTRF Prostaglandin E2 kit - 500 pts

Product Number : 62P2APEG Lot Number : 14D

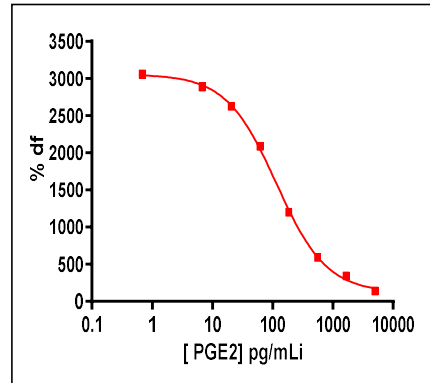
Expiration date : 2025-04-05 Document version : 1

Reagents :	Lot n°
HTRF Diluent 2 - 20 mL	10RA
HTRF Detect. Buf. 2 - 7 mL	17RA
HTRF Prostaglandin E2 kit - 500 pts d2 reagent	14RA
HTRF Prostaglandin E2 kit - Std	12RA
HTRF Prostaglandin E2 kit - 500 pts Eu Cryptate antibody	14RA

Results :

Incubation : Overnight at 4°C
 Plate : 384 wells small volume white plate 20 µL
 Reader : Pherastar Plus
 Fitting model : nonlinear regression, sigmoidal dose response (variable slope)

PGE2	Conc. (pg/ml)	DF %
Std 0	0	3057
Std 1	6,85	2887
Std 2	20,5	2623
Std 3	61,7	2087
Std 4	185,1	1196
Std 5	555,5	592
Std 6	1666,6	340
Std 7	5000	138



HTRF Prostaglandin E2 kit - 500 pts - Lot n° : 14D

Accepted Rejected

Quality control date : 2023-12-08	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.