

HTRF® Research Reagents

HTRF Prostaglandin E2 kit - 10K pts

Product Number : 62P2APEH Lot Number : 10X

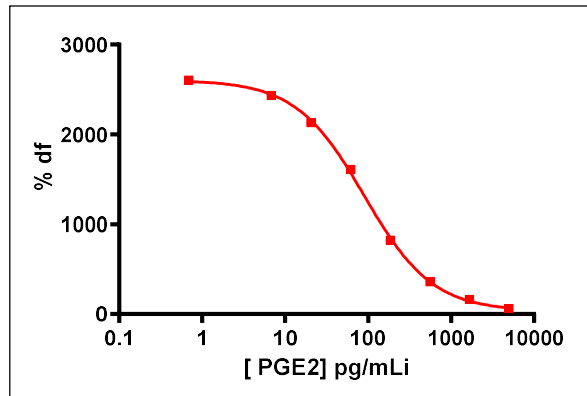
Expiration date : 2025-11-30 Document version : 1

Reagents :	Lot n°
HTRF Diluent 2 - 20 mL	12A
HTRF Detect. Buf. 3 - 105 mL	29RA
HTRF Prostaglandin E2 kit - 10K pts d2 reagent	09RA
HTRF Prostaglandin E2 kit - Std	13A
HTRF Prostaglandin E2 kit - 10K pts Eu Cryptate antibody	10RA

Results :

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

PGE2	Conc. (pg/ml)	DF %
Std 0	0	2605%
Std 1	6,85	2434%
Std 2	20,5	2130%
Std 3	61,7	1611%
Std 4	185,1	821%
Std 5	555,5	362%
Std 6	1666,6	163%
Std 7	5000	60%



Assay characteristics :

	Found value	Specifications
Delta F%	2605%	2600 - 3700 %
Detection limite	3,9 pg/ml	≤ 15 pg/ml
EC 50	90	90-140 pg/ml

HTRF Prostaglandin E2 kit - 10K pts - Lot n° : 10X

Accepted

Rejected

Quality control date : 2024-10-02	Supervisor's initials : MP
-----------------------------------	----------------------------

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.