

HTRF® Research Reagents

HTRF LTC4 kit - 500 pts

Product Number : 64LC4PEG Lot Number : 14E

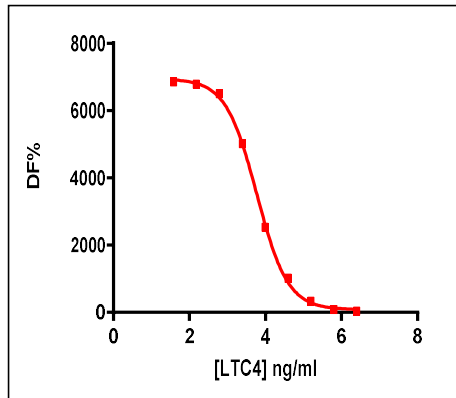
Expiration date : Document version : 1

Reagents :	Lot n°
HTRF Diluent 3 - 20 mL	13AR
HTRF Detect. Buf. 3 - 7 mL	30AR
HTRF LTC4 kit - Std	09RA
HTRF LTC4 kit - 500 pts d2 reagent	14RC
HTRF LTC4 kit - 500 pts Eu Cryptate antibody	14RB

Results :

Incubation : 1h at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : Pherastar Plus
 Fitting model :

LTC4	Conc. (pg/ml) _i	DF (%)
Std 0	0	6848%
Std 1	0,038	6785%
Std 2	0,152	6501%
Std 3	0,61	5018%
Std 4	2,44	2533%
Std 5	9,76	1012%
Std 6	39,06	326%
Std 7	156,25	85%
Std 8	625	29%
Std 9	2500	15%



Assay characteristics :

	Found value	Specifications Rev 01
DF% max	6848%	5900% - 8900%
EC50	1,5 ng/ml	1.5-3.2 ng/ml

HTRF LTC4 kit - 500 pts - Lot n° : 14E

Accepted

Rejected

Quality control date : 2024-02-19	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.