

HTRF® Research Reagents

HTRF CHO HCP Kit - 10K pts

Product Number : 64CHOPEH Lot Number : 02A

Expiration date : 2025-06-12 Document version : 1

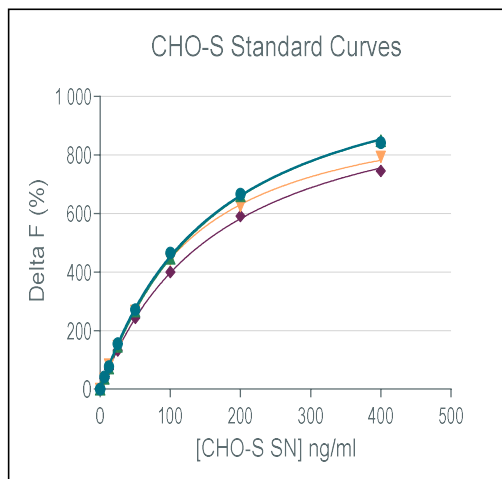
Reagents :	Lot n°
HTRF Diluent 5 (5X) - 10 mL	21RA
HTRF Detect. Buf. 17 - 50 mL	01RA
HTRF CHO HCP Kit - 10K pts d2 antibody	02A
HTRF CHO HCP Kit - 10K pts Eu Cryptate antibody	02A
HTRF CHO HCP Kit - Std	01RB

Results :

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

Assay characteristics :

CHO	Conc (ng/ml)	Delta F %	Delta F% Accepted Range Provisoire
Std 0	0	0%	0-0
Std 1	6,25	37%	25 -37
Std 2	12,5	69%	51 - 77
Std 3	25	133%	99 - 149
Std 4	50	244%	182 - 273
Std 5	100	400%	314 - 470
Std 6	200	591%	484 - 727
Std 7	400	745%	630 - 946



HTRF CHO HCP Kit - 10K pts - Lot n° : 02A

Accepted

Rejected

Quality control date : 2024-02-09 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