

HTRF® Research Reagents

HTRF Methyltransferase kit - 10K pts

Product Number : 62SAHPEH Lot Number : 12M

Expiration date : 23/06/2024 Document version : 1

Reagents :	Lot n°
HTRF Methyltransferase kit - Std	12RA
HTRF Epig. diluent buf. - 13 mL	10RA
HTRF Methyltransferase kit - 10K pts modop	3
HTRF Methyltransferase kit - 10K pts d2	12RA
HTRF Methyltransferase kit - 10K pts Tb Cryptate antibody	12RA
HTRF Epig. Detect. Buf. 1 - 24 mL	15RA
HTRF Epig. Detect. Buf. 2 - 130 mL	13RA
HTRF Ultrapure SAM cofactor - 2.4 µmol	16RA

Results :

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

Assay characteristics :

Norm 20µM	Found value	Specifications
IC50	45,2	30 - 60 nM
S/B	13,3	12 18

Norm 0.5 µM	Found value	Specifications
IC50	8,2	7 - 14 nM
S/B	9,5	6 12

HTRF Methyltransferase kit - 10K pts - Lot n° : 12M

 Accepted

 Rejected

Quality control date : 2022-03-24	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.