

HTRF® Research Reagents

HTRF Human SMN kit - 500 pts

Product Number : 63ADK033PEG Lot Number : 14C

Expiration date : 2026-11-29 Document version : 1

Reagents :	Lot n°
HTRF Detect. Buf. 3 - 7 mL	30AR
HTRF Human SMN kit - 500 pts d2 antibody	14A
HTRF Human SMN kit - 500 pts Eu Cryptate antibody	14A
HTRF Human SMN kit - Ctrl lysate	14A
HTRF P-T prot. - Lysis Buf.2 (4X) 2 mL	13AR

Results :

Incubation : Overnight at RT
 Plate : 384 wells small volume white plate 20 µL
 Reader : /
 Fitting model : /

Assay characteristics :

	Found value	Specifications Rev 01
S/B	4.34	>4

HTRF Human SMN kit - 500 pts - Lot n° : 14C

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