

HTRF® Research Reagents

HTRF AAV1 Capsid Kit - Std
Product number: 64AAV1CDA

Lot number: 01A

Product characteristics:

- **Packaging Unit** 3 tests/vial
- **Conditioning buffer** PBS + 0.014% Tween
- **Appearance** Frozen form
- **Storage conditions** -80 °C
- **Expiration date** 2026.03.15
- **Volume per vial** 90.00 µL
- **Active Moiety Concentration (St)** 4.99999998E11 VP/ml

Weight indications:

	Quantity per vial	MW (Da)
Active Moiety (St)	4.999999898E10 VP	N/A

Please note

Conc St en VP/mL et non mg/ml

Qt St par tube en VP et non mg

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.