

AAV Reference Standards

rAAV8 full capsids – Reference standard material

Product number: IP24016 Lot Number: AAV8-10251_V1

Material provided: rAAV8-CMV-EGFP-WPRE

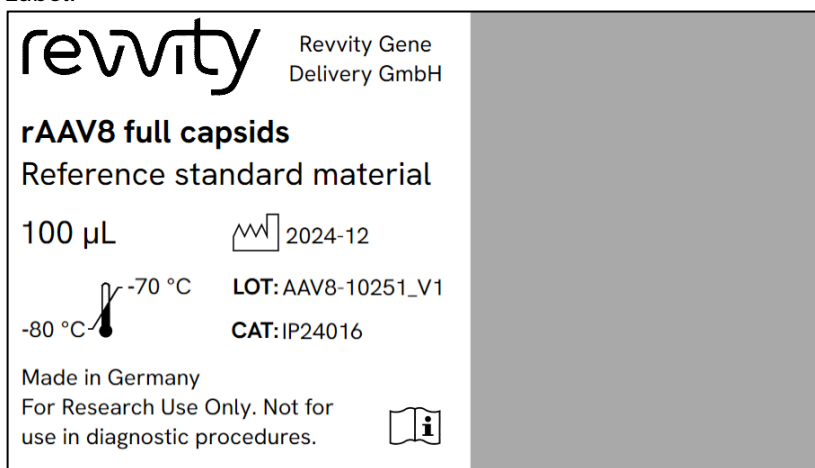
Format: 100 µL

Manufacturing date: 12/17/2024 Document version: 1

Product characteristics:

- Packaging unit Transparent polypropylene screw-cap tube with EPDM rubber O-ring
- Conditioning buffer DPBS + 0.001% Poloxamer 188
- Appearance Frozen liquid
- Storage conditions Store from -70 °C to -80 °C, do not aliquot, avoid freeze-thaw cycles
- Stability Tested yearly
- Transgene CMV-EGFP-WPRE flanked by AAV2 ITRs
- Biosafety level BSL 1
- Particle concentration >1E12 particles/mL
- Endotoxin <1 EU/mL
- Mycoplasma Negative
- Theoretical molecular weight Approximately 4.9 MDa

Label:



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.