

Protocol • Lenti-Pac™ Lentivirus Storage Buffer • Cat. No. LT039

Description

GeneCopoeia has developed **Lenti-Pac™ Lentivirus Storage Buffer**, which can improve the stability of the lentivirus. Using this product to dissolve lentivirus particles can improve the stability of lentivirus at -80° C and effectively preserve functional viral particles during repetitive freeze-thaw cycles, effectively slowing down the decline in virus titer and preserving lentiviral infectivity. It is suitable for resuspending lentivirus after ultracentrifugation, polyethylene glycol precipitation (PEG precipitation), and ultrafiltration concentration.

- Significantly improve the freeze-thaw stability of lentivirus.
- Effectively preserve lentiviral infectivity.
- Performance.
- Suitable for stable storage of any lentiviral particles.

Content and Storage

Component	Volume	Shipping Condition	Storage Condition
Lenti-Pac™ Lentivirus Storage Buffer	20 ml	Room temperature	2~8°C Stable for at least 12 months

Protocol

Resuspension of lentivirus after ultracentrifugation, PEG precipitation, and ultrafiltration concentration:

1. According to the expected concentration multiple, add an appropriate amount of lentivirus storage buffer (for example, if 10 mL of lentivirus stock solution is expected to be concentrated 100 times, add 100 µl of lentivirus storage buffer), and then use a pipette to gently blow up the virus precipitate.
2. Aliquot lentivirus and store at -80°C.