

MANUAL

LentiBOOST™

Transduction Enhancer

Part number: LentiBOOST™-P

Storage: -15°C to -25°C

Shipped at room temperature

Version: 1

Date: July 2024

THE USE OF THE PRODUCT IS SUBJECT TO THE LIMITED USE LABEL LICENSE AGREEMENT AS SET OUT AT THE END OF THIS MANUAL.

1. PRODUCT DESCRIPTION

LentiBOOST™ technology enhances the uptake of lentiviral vectors into mammalian cells. Compared to commonly used transduction enhancers like polybrene or protamine sulfate, it does not negatively affect cell viability or cell growth.

Furthermore, for human CD34+ hematopoietic stem cells (section 6) and progenitor cells studies have indicated that LentiBOOST transduction enhancer (TE) allows to maintain the differentiation potential of the cells after lentiviral transduction (Hauber et al. 2018). The authors observed high transduction rates while limiting the vector copy number (VCN close to 5) to a moderate increase that reaches each individual cell as requested by regulatory guidelines.

For human T-cells (section 7) a group of scientists has shown that the use of LentiBOOST TE technology for lentiviral transduction can result in a 2-fold increase in transduction rate, without observing a cytotoxic effect on the cells (Simon et al. 2019).

LentiBOOST technology is therefore an excellent tool for enhancing lentiviral transduction of sensitive primary cells including hematopoietic stem cells and T-cells.

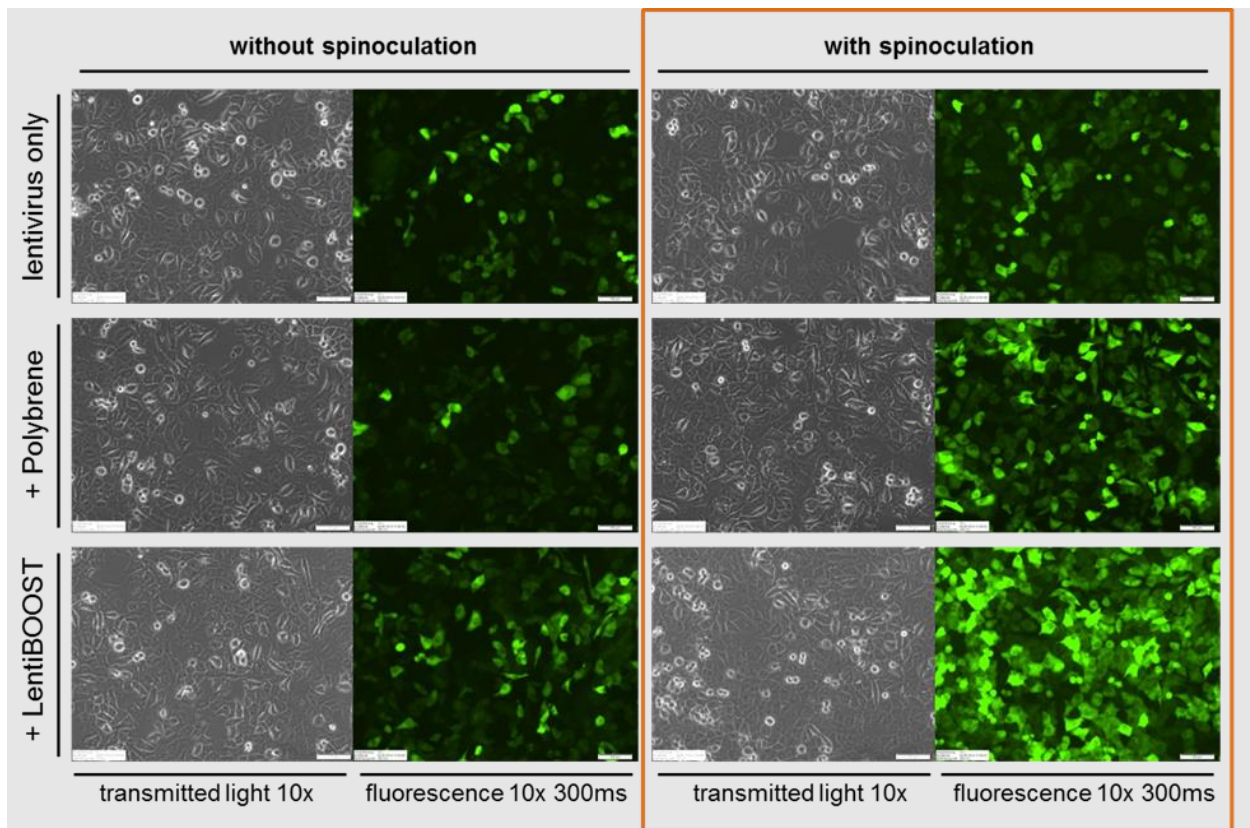


Fig. 1: Lentiviral transduction efficiencies in H1299 48h after transduction with lentivirus LV-CMV-GFP. Protocols according to the instructions of this manual.

2. LentiBOOST TE pharma grade non GMP

LentiBOOST-P TE 100 mg/ml aqueous solution

3. Expiration Date

Not to be used after expiration date.

4. Storage

Store at -15°C to -25°C.

5. Thawing and handling

It is recommended to thaw LentiBOOST enhancer between +4°C and +25°C.

Before opening the tube, spin down briefly to remove any liquid from the lid. To keep for later use, you can aliquot LentiBOOST enhancer using aseptic technique and store at -15°C to -25°C. LentiBOOST TE can be stored at +4°C for 1 month. No loss of functionality* was observed after 5 freeze-thaw cycles.

*based on an internal standardized assay, no conclusion can be drawn to any customer specific cell types.

PLEASE NOTE: The protocols given below are suggestions only based on (published) customer data. Protocols must be adapted depending on customer-specific conditions.

6. Transduction Protocol for CD34+ Transduction

Preparation of human CD34+ HSC

CD34+ hematopoietic stem cells (HSC) are isolated from PBMC using standard protocols.

Day 1: Seeding cells

CD34+ HSC are cultured according to standard protocols. Cells are incubated at 37°C in a humidified incubator containing 5% CO₂.

Day 2: Transduction

The following day, CD34+ HSCs are counted and 1×10^6 cells are seeded per well in a 24 well plate. LentiBOOST enhancer and lentiviral vector are directly added to the cells.

For an initial experiment, it is recommended to use MOIs between 2 - 30 for transduction and to add LentiBOOST TE at an initial concentration of 1 mg/ml (1:100) of the total volume (medium + virus). In a second experiment, it is recommended to titrate LentiBOOST enhancer in the range of 5 mg/ml – 0.1 mg/ml (1:20-1:1000) to determine the minimal active concentration.

Optionally spinoculation can be performed at 600g for 90 minutes at room temperature.

Day 3: Medium exchange

- Aspirate medium from transduced cells and add appropriate amount of normal growth medium.

7. Transduction Protocol T-cells

Seeding and pre-stimulation of cells

- Primary T Cells are stimulated according to standard protocols. This step should be adapted according to individual protocols.

Day 1: Transduction

- Thaw lentivirus at +4°C.
- Prepare 500µl of culture medium and add LentiBOOST enhancer at 1 mg/ml (1:100) as starting point.
- Concentrations of LentiBOOST TE can be tested between 0.1 mg/ml and 5 mg/ml.
- Add viral vector to medium at desired MOI and mix gently.
- For an initial experiment, it is recommended to use MOIs between 2-30.
- Pellet 10^6 cells by centrifugation (cell number can be adapted according to needs).
- Mix the cell pellet with the prepared medium containing LentiBOOST TE and viral vector.
- Seed the cells into a 24 well plate.
- Spinoculation (optional): Centrifuge cell culture plate for 90 min at 800 g at RT.
- Incubate the overnight at 37°C and 5% CO₂.

Day 2: Medium change

- Exchange medium according to standard protocols.

8. Literature

Hauber, I., Beschorner, N., Schrödel, S., Chemnitz, J., Kröger, N., Hauber, J., & Thirion, C. (2018). Improving Lentiviral Transduction of CD34+ Hematopoietic Stem and Progenitor Cells. Human Gene Therapy Methods, hgtb.2017.085. <https://doi.org/10.1089/hgtb.2017.085>

Simon, B., Harrer, D. C., Thirion, C., Schuler-Thurner, B., Schuler, G., & Uslu, U. (2019). Enhancing lentiviral transduction to generate melanoma-specific human T cells for cancer immunotherapy. Journal of Immunological Methods, 472(April), 55–64. <https://doi.org/10.1016/j.jim.2019.06.015>

9. Troubleshooting

Here are some general approaches that LentiBOOST TE users have found useful when troubleshooting issues. However, please note that users will need to account for the specific system.

Trouble	Possible reason	Solution
Low transduction efficiency	MOI used was too low	Use higher amounts of lentivirus up to MOI 50
	Cells are very hard to transduce	Include spinoculation step 800 g for up to 90 min at room temperature (in cell culture plates)
		Increase concentration of LentiBOOST enhancer up to 5 mg/ml (1:20)
		Add Protamine sulfate at 5 µg/ml additionally to LentiBOOST TE
Low viability	Cells are sensitive to LentiBOOST enhancer	Decrease concentration of LentiBOOST TE to e.g. 0.2 mg/ml, 0.1 mg/ml (1:500, 1:1000)
	Cells are sensitive to spinoculation	Try protocol without spinoculation Reduce duration Reduce velocity
	Cells are sensitive to lentiviral vectors	Change medium 4 h after transduction or directly after centrifugation

Limited Use Label License Agreement

BEFORE USING THIS PRODUCT, PLEASE READ THE TERMS AND CONDITIONS SET FORTH IN THIS LIMITED USE LABEL LICENSE AGREEMENT (THE "AGREEMENT"). YOUR USE OF THIS PRODUCT SHALL CONSTITUTE ACKNOWLEDGMENT AND ACCEPTANCE OF THESE TERMS AND CONDITIONS.

If you are not willing to use the Product pursuant to the terms and conditions of this ("Agreement"), please contact Revvity Gene Delivery GmbH to return the unused and unopened Product for a full credit. Address: Revvity Gene Delivery GmbH, Head of Logistics, Am Haag 6, 82166 Gräfelfing, Germany, Email: mun.info@revvity.com, Tel.: +49 89 70096199-0

IMPORTANT INSTRUCTIONS – READ CAREFULLY: This Agreement is the legally binding agreement between you (hereinafter "**Licensee**") and Revvity Gene Delivery GmbH ("**Revvity Gene Delivery**") (individually, a "**Party**" or together, the "**Parties**"), for use of LentiBOOST™ TE pharma grade non GMP ("**Product**"), whether such Product is provided by Revvity Gene Delivery, its affiliates, (sub)licensees, resellers, or distributors. If the use of the Product is subject to multiple limited use label licenses, the terms of the most restrictive limited use label license shall control. **Licensee's use of the Product shall constitute acknowledgment and acceptance of the terms and conditions of this Agreement.**

1. Use of Products by Licensee. Revvity Gene Delivery grants to Licensee a non-exclusive, non-transferable, non-sublicensable (other than in compliance with Section 5 below), limited right to use only the purchased amount of the Product for Research Use Only, subject, and provided that (i) no further license, right or permission is granted hereunder, and (ii) not for the Excluded Uses (as defined below). **"Research Use Only"** means internal laboratory research by Licensee for the sole benefit of Licensee, and specifically excluding use (a) for therapeutic, diagnostic, vaccine or prophylactic purposes, and (b) any applications which require regulatory approval, including any in vitro diagnostic or therapeutic applications in humans, any ex vivo use to enable gene therapy, or (c) for any in vivo use in humans for any purpose.

2. Excluded Uses. The Product is not licensed, and Licensee may not use, the Product for any Commercial Purpose, Development Purpose or in the Excluded Field, and no licenses are granted expressly, by implication, by estoppel, or otherwise for any such Commercial Purpose, Development Purpose or in the Excluded Field. "**Commercial Purpose**" means any commercial activity for a third party for consideration, including but not limited to: (a) any use, directly or indirectly, in manufacturing and/or production, or quality control; (b) any use to provide a service, information, or data; (c) any sale, resale, leasing, distribution, transfer, licensing, importation of, or other transfer of the Product (whole or in part) or derivatives thereof to any third party, in all cases above whether or not such transfer is limited for Research Use Only. "**Development Purpose**" means any clinical activity following IND-enabling preclinical toxicological studies or equivalents thereof. "**Excluded Field**" means any use of the Product for ex vivo gene therapy for hemoglobinopathies. Licensee acknowledges and agrees that Licensee shall not have the right to authorize any third party to use or sell the Product or derivatives of the Product which contain poloxamers protected (other than in compliance with Section 5 below) by the below patent

applications/patents and other intellectual property rights owned or controlled by Revvity Gene Delivery or its affiliates.

3. Licensor's Patent Protection and Rights to Inventor's Inventions. The Product and its use are protected under patent applications/patents (WO2013127964) and other intellectual property rights owned or controlled by Revvity Gene Delivery or its affiliates. Revvity Gene Delivery and its affiliates retain all rights not expressly granted herein, and there are no implied licenses or rights granted herein by estoppel, implication or otherwise.

If Licensee's use of the Product as permitted under this Agreement leads to any invention by Licensee that does not incorporate the Product and does not require the use of the Product, then as between Revvity Gene Delivery and Licensee, Licensee shall own all such inventions other than Improvements and shall have the first right to file any patent application covering such inventions at its own cost and expense. **"Improvement"** means any further development, enhancement or improvement by either Party that is within the scope of the disclosure contents of the published PCT application WO 2013/127964 A1, and, without limitation the generality of the foregoing, (i) modifications of concentrations of substances, compounds such as poloxamers and secondary agents such as polycationic polymers and polycationic peptides, (ii) combinations of compounds or agents disclosed in the above application with further agents, or (iii) uses of compounds or agents disclosed in the above application for medical indications; and such Improvement shall be, remain, and become Revvity Gene Delivery's sole and absolute property.

For clarity, patent applications permitted by this Section 3 may describe the Product but shall not include any claim (i) reciting the Products or use thereof, (ii) covering the Products or use thereof, or (iii) otherwise excluding others from making, using, or selling the Products.

In the event Licensee files in any jurisdiction any application for a patent having any claim prohibited under this Section 3, in addition to any other remedies available to Revvity Gene Delivery, Licensee hereby automatically grants to Revvity Gene Delivery for so long as one or more claims prohibited under this Section 3 remain in such application or patent, a non-exclusive, irrevocable, fully paid-up license (with the right to sub-license also in multiple tiers) to research, have researched, develop, have developed, manufacture, have manufactured, use, have used, import, have imported, offer to sell, have sold, commercialize and/or have commercialized subject matter claimed in any and all patents that issue. Such license to Revvity Gene Delivery shall not include the right to use Licensee's intellectual property that is separable or independent of the Product, including but not limited to Licensee's background intellectual property.

4. Protection of Trade Secrets. Licensee acknowledges and agrees that the Product may contain valuable trade secret information and other proprietary information. Accordingly, the Parties shall not, at any time during the term of this Agreement or thereafter, reverse engineer or

otherwise attempt to discern the trade secret information or other proprietary information, nor shall the Parties permit any Third Party to do any of the foregoing.

5. Use by Licensee's Third-Party Contractors. Licensee shall be permitted to provide the Product to a third party contractor to use the Product for the sole benefit of the Licensee and its internal laboratory research, provided however that a legal agreement is in place and in force between Licensee and such third party contractor, whose terms will provide that the ownership of the Product is not transferred to such third party contractor and Licensee can require the return of the Product at any time.

6. Compliance. Licensee shall use (and cause its third party contractors to use) the Product at all times in compliance with all applicable laws and regulations, including applicable human health and animal welfare laws and regulations. Any violation of this stipulation shall constitute a breach of this Agreement.

7. Misuse of the Products and Indemnification terms. Licensee will defend, indemnify and hold harmless Revvity Gene Delivery and its affiliates, and its and their managers, directors, officers, employees, sponsors, and agents (collectively the "**Indemnified Parties**") against any and all (alleged) (i) liability, (ii) loss, (iii) damage, (iv) costs, (v) claims, or (vi) expenses, including (patent) attorneys' fees (collectively the "**Indemnified Losses**"), all arising out of or in connection with this Agreement, including, without limitation Indemnified Losses resulting from any use of the Product (whole or in part) and any derivatives thereof by the Licensee, Licensee's employees, representatives, students, consultants, third party contractors or agents. Licensee will indemnify and hold harmless the Indemnified Parties against any and all Indemnified Losses resulting from, arising out of or relating to: (i) product liability claims of any nature; (ii) claims arising from Licensee's failure to comply with governmental regulations relating in any way to use or storage of the Product (whole or in part) and any derivatives thereof; (iii) Licensee's breach of this Agreement; and (iv) claims by a third party that Licensee, Licensee's representatives, employees', representatives', students', third party contractors' or agents' use of the Product infringes or violates any patent, copyright, trademark or property rights of such third party.

8. **DISCLAIMER OF WARRANTY.** NEITHER REVVITY GENE DELIVERY NOR ANY OF ITS AFFILIATES MAKES ANY WARRANTY, EXPRESS, OR IMPLIED WITH RESPECT TO THE PRODUCT, INCLUDING ANY WARRANTY OF MERCHANTABILITY OR FITNESS FOR ANY PARTICULAR PURPOSE OR THAT THE PRODUCT DOES NOT INFRINGE ANY PATENT, COPYRIGHT, TRADEMARK, OR OTHER RIGHTS OR ANY OTHER EXPRESS OR IMPLIED WARRANTIES.

9. Term. This Agreement shall commence upon the receipt of the Product and, subject to Section 10, continue as long as the Product is in the possession of the Licensee.

10. Termination. Either Party may terminate this Agreement for any reason upon thirty (30) days' prior written notice to the other Party. The rights and obligations under Sections 3, 4, 6, 7, and 8 shall survive any termination, expiration, or completion of this Agreement with respect to information generated and activities and events occurring prior thereto. Upon expiration or any

termination of this Agreement, Licensee shall promptly destroy all remaining Product, any derivatives thereof, and all material made using the Product.

11. Entire Agreement and Assignability. This Agreement sets forth the complete and entire agreement of the Parties with respect to the Product and supersedes and terminates all prior agreements and understandings between the Parties with respect to the Product. No subsequent amendment or addition to this Agreement shall be binding upon the Parties unless reduced to writing and signed by the respective authorized representatives of the Parties. This Agreement shall not be assigned or otherwise transferred by Licensee.

12. Governing Law and Jurisdiction. This Agreement shall be governed by and construed in accordance with the laws of Germany, excluding its conflict of laws principles. For all disputes arising out of or in connection with this Agreement or its validity the Parties submit to the exclusive jurisdiction of the courts having jurisdiction for the seat of Licensor.

Additional Obligations

Licensee will promptly inform Revvity Gene Delivery in writing at the address specified below in the event of any uses in conflict with those permitted by this Agreement.

For information on purchasing a license to the Product for purposes other than Licensed Research Use, please contact Revvity Gene Delivery:

Business Development & Licensing

Revvity Gene Delivery GmbH

Street: Am Haag 6

Postal code: 82166

City: Gräfelfing

Country: Germany

Email: mun.licensing@revvity.com

Tel.: +49 89 70096199-0

Revvity Gene Delivery reserves the right to grant, or not to grant, permission for any such other uses.



The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application.

www.revvity.com

revvity

Revvity Gene Delivery GmbH
Am Haag 6
82166 Gräfelfing, Germany

(800) 762-4000
www.revvity.com

For a complete listing of our global offices, visit www.revvity.com
Copyright ©2025, Revvity, Inc. All rights reserved.