

Nupherin™

Catalog #: BML-SE225

ACTIVITY: The Nupherin™ reagent is a specially designed peptide formulation that serves as an enhancer of plasmid transfer and expression in lipofected cell lines. The Nupherin™ reagent is used with cationic lipid and is compatible with most commercially available cationic lipid formulations.

Use at 5 to 50µg Nupherin™ reagent per 1µg of plasmid.

STORAGE: Nupherin™ is supplied in sterile DNase-free distilled water, frozen at 3 mg/ml.

Store frozen at -80°C.

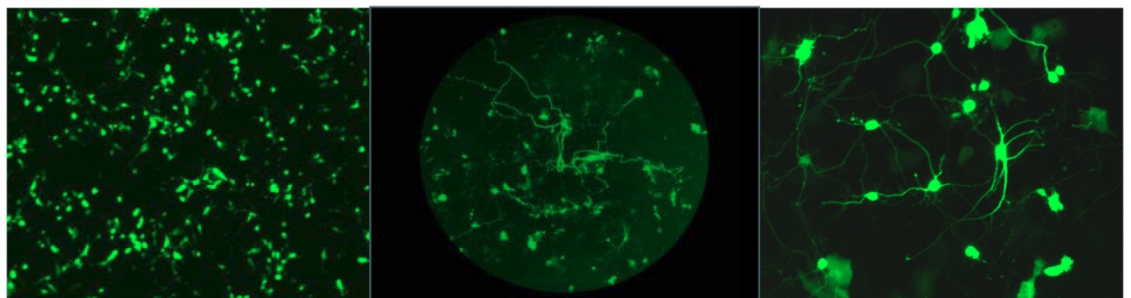
On first use, aliquot the reagent and store at -80°C.

Thawed aliquots of Nupherin™ can be refrozen after use for long term storage or refrigerated for up to 1 week. Do not dilute the reagent.

Avoid repeated freeze/thaw cycles.

REFERENCES:

1. A. Subramanian, et al. Nature Biotechnology 1999 17 873
2. N.B. Arnold, et al. Cancer Res. 2004 64 3599
3. L.J. Karai, et al. J. Biol. Chem. 2004 279 16377
4. V. Chennathukuzhi, et al. Proc. Natl Acad Sci (USA) 2003 100 15566
5. E. Trushina, et al. Proc. Natl Acad Sci (USA) 2003 100 12171
6. A. Devillers-Thiery et al. Biol Cell 2003 95 373
7. R. Nashmi, et al. J Neurosci. 2003 23 11554
8. E.M. Slimko, et al. J Neurosci Methods 2003 124 75
9. E.M. Slimko, et al. J. Neurosci. 2002 22 7373



A.

B.

C.

Figure A: Chicken Embryonic Retinal Ganglia

Figure B: Primary neuron lipofection

Figure C: Embryonic rat primary hippocampal neuron

**FOR RESEARCH USE ONLY.
NOT FOR USE IN DIAGNOSTIC PROCEDURES.**

SAMPLE PROTOCOL FOR NUPHERIN™

1. Culture cells on glass or plastic coverslips placed in 6-well (9.6cm²) or 24-well (2.0cm²) plates or culture cells directly in 6-well or 24-well plates. Cells and cell lines can be transfected at subconfluence, confluence, or in a post-mitotic state.
2. TUBE A: For each well of a 24-well plate to be transfected, use 0.5 to 3µg plasmid DNA per well. 1µg or less of DNA is preferred and may result in higher transfection rates. Dilute plasmid into 150µl of serum-free phenol red-free DMEM or OPTIMEM™ in Tube A. Add 3 to 20µl of the Nupherin™ reagent (9 to 60µg) per µg plasmid to Tube A. Gently mix by pipetting (do not vortex). Incubate 15 min at room temperature before mixing with Tube B.
3. TUBE B: Add cationic lipid (e.g. 1 to 4µl Lipofectamine™ at 2µg/µl) to a final volume of 150µl serum-free phenol red-free DMEM or OPTIMEM™ in Tube B. For other cationic lipid reagents, use: 0.5-3µl Fugene™, 1-3µl lipofectin™, 1-4µl lipofectamine™, 2-6µl Geneporter™.
4. Gently mix Tube A and Tube B by pipetting (do not vortex). Incubate for 40 minutes at room temperature.
5. Rinse cells in serum-free medium, overlay cells in each 2-cm² well with the transfection solution (about 300µl).
6. Centrifuge the plate at 100g (1000 rpm in swinging bucket rotor) for 3 to 5 min at room temperature.
7. Incubate the plate at 37°C for 0.5 to 4 hr. Replace with serum-containing growth medium.

APPLICATION EXAMPLE

Recommended Protocol for Post-mitotic Primary Neurons

| | 2.0cm² well (24-well plate) | 9.6cm² well (6-well plate) |
|----------------------------|---|--|
| plasmid (1µg/µl) | 1µl | 2µl |
| Nupherin (3µg/µl) | 10µl | 20µl |
| OPTIMEM™(1X) | <u>140µl</u> 150µl Tube A | <u>380µl</u> 400µl Tube A |
| Lipofectamine™ (2µg/µl) | 1µl | 3µl |
| OPTIMEM™ (1X) | <u>150µl</u> 150µl Tube B | <u>400µl</u> 400µl Tube B |

Gently overlay 300µl of transfection mixture (Tube A + Tube B) over 2.0cm²-well or 800µl over 9.6 cm²-well, centrifuge 100g for 3 min, incubate 30 min, and replace with growth media.

All trademarks are the property of their respective owners. Nupherin™ is a trademark of Promid, LLC. Lipofectamine™, OPTIMEM™ are trademarks of Life Technology, Inc. Nothing in this document should be construed as an authorization or implicit license to make or sell any components covered under any patent. Use of this product for clinical diagnostic, human, or veterinary applications is strictly forbidden.

USE FOR RESEARCH PURPOSES ONLY

Unless otherwise specified expressly on the packaging, all products sold hereunder are intended for and may be used for research purposes only and may not be used for food, drug, cosmetic or household use or for the diagnosis or treatment of human beings. Purchase does not include any right or license to use, develop or otherwise exploit these products commercially. Any commercial use, development or exploitation of these products or development using these products without the express written authorization of Enzo Life Sciences, Inc. is strictly prohibited. Buyer assumes all risk and liability for the use and/or results obtained by the use of the products covered by this invoice whether used singularly or in combination with other products.

LIMITED WARRANTY; DISCLAIMER OF WARRANTIES

These products are offered under a limited warranty. The products are guaranteed to meet all appropriate specifications described in the package insert at the time of shipment. Enzo Life Sciences' sole obligation is to replace the product to the extent of the purchasing price. All claims must be made to Enzo Life Sciences, Inc., within five (5) days of receipt of order. THIS WARRANTY IS EXPRESSLY IN LIEU OF ANY OTHER WARRANTIES OR LIABILITIES, EXPRESS OR IMPLIED, INCLUDING WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, AND NON- INFRINGEMENT OF THE PATENT OR OTHER INTELLECTUAL PROPERTY RIGHTS OF OTHERS, AND ALL SUCH WARRANTIES (AND ANY OTHER WARRANTIES IMPLIED BY LAW) ARE EXPRESSLY DISCLAIMED.

TRADEMARKS AND PATENTS

Several Enzo Life Sciences products and product applications are covered by US and foreign patents and patents pending. Enzo is a trademark of Enzo Life Sciences, Inc.

GLOBAL HEADQUARTERS

Enzo Life Sciences Inc.
10 Executive Boulevard
Farmingdale, NY 11735
Toll-Free: 1.800.942.0430
Phone: 631.694.7070
Fax: 631.694.7501
info-usa@enzolifesciences.com

EUROPE/ASIA

Enzo Life Sciences (ELS) AG
Industriestrasse 17
CH-4415 Lausen
Switzerland
Phone: +41/0 61 926 89 89
Fax: +41/0 61 926 89 79
info-ch@enzolifesciences.com

For local distributors and detailed product information visit us online:
www.enzolifesciences.com