

Nupherin™

Transfection Reagent

Greatly enhances transfection efficiency

Nupherin™ dramatically enhances the performance of commercially available cationic lipids to allow high-efficiency lipofection (10-40%) of neuronal, ganglion and other difficult to transfect cells. Efficient lipofection is limited by the ability of transfected DNA to enter the nucleus – a particular obstacle in nondividing or slow dividing cells like neurons. Nupherin, a peptide based reagent that combines a non-classical nuclear localization sequence (M9) with a nucleotide binding domain, helps to transport DNA into the transfected cell's nucleus. Nupherin has been demonstrated to provide high efficiency lipofection of primary rat hippocampal neurons, chicken embryonic retinal ganglionic cells, rat pheochromocytoma (PC12) cells and other recalcitrant cells. Fully compatible with cells grown on glass coverslips, in plastic plates, or in tissue-culture flasks.

Citations: 22

[View Online »](#)

Ordering Information

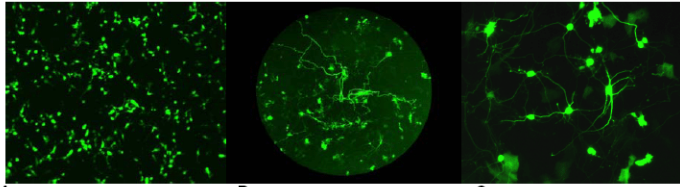
[Order Online »](#)

BML-SE225-0075	0.75mg
BML-SE225-0003	3mg

Manuals, SDS & CofA

[View Online »](#)

- Dramatically enhances performance of commercially available cationic lipids
- Allows high-efficiency lipofection (10-40%) of neuronal and ganglion cells
- Fully compatible with cells grown on glass coverslips, plastic plates or tissue-culture flasks



A.

B.

C.

Figures: **A.** Chicken Embryonic Retinal Ganglia, **B.** Primary neuron lipofection, **C.** Embryonic rat primary hippocampal neuron

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. On first use, aliquot the reagent and store at -80°C. Thawed aliquots of Nupherin™ can be frozen after use for long term storage or refrigerated for up to 1 week. Do not dilute reagent.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Formulation	Liquid. Sterile solution.
Quantity	The 3 mg size provides enough reagent for 100-200 transfections.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)