

LyoVec™

A cationic lipid for transfecting mammalian cells

Catalog code: lyec-1, lyec-2, lyec-3

<https://www.invivogen.com/transfection-reagents>

For research use only

Version 21121-MM

PRODUCT INFORMATION

Contents:

LyoVec™ is supplied as a lyophilized lipid formulation to be reconstituted with water. It is compatible with all cell culture media. This product is available in 3 pack sizes:

- **lyec-1:** 5 small vials (125 µg per vial), sufficient to perform 200 reactions
- **lyec-2:** 10 small vials (125 µg per vial), sufficient to perform 400 reactions
- **lyec-3:** 2 large vials (500 µg per vial), sufficient to perform 320 reactions

One reaction is a transfection of 2-4 x 10⁵ cells/well in a 12-well plate.

Storage and stability:

- LyoVec™ is shipped at room temperature and should be stored at 4 °C. Lyophilized product is stable for 1 year when properly stored.
- After resuspension, store at 4 °C. Resuspended LyoVec™ is stable for 6 months when properly stored.

Quality control:

- Transfection efficiency of LyoVec™ has been confirmed in HEK293-derived cells.
- Cytotoxicity has been assessed in HEK293-derived cells.

DESCRIPTION

LyoVec™ is a proprietary cationic lipid belonging to the phosphonolipid family, originally described by Floch *et al.*, as an efficient transfection reagent both *in vitro* and *in vivo*^{1,2}. The major constituent of LyoVec™ is the phosphonolipid DTCPTA, which is coupled to DiPPE, a neutral lipid that helps destabilize membrane bilayers and increases the *in vitro* transfection efficiency of LyoVec™. The positive charge of the phosphonolipid enables LyoVec™ to bind to plasmid DNA, where the structure promotes fusion with cellular membranes for efficient DNA delivery. LyoVec™ is designed as a transfection reagent for mammalian cells.

Additionally, LyoVec™ can be used as a nucleic acid complexing agent to facilitate the cellular entry of RNA or DNA-based oligonucleotides, such as RIG-I ligands (e.g. Poly(dA:dT)³ and poly(I:C)⁴; see Related Products on the next page) and TLR7/8 ligands (e.g. ssPolyU⁵). This complexation step is crucial for inducing an effective response to the nucleic acids by intracellular pathogen recognition receptors.

1. Flochet *et al.*, 1997. Cationic phosphonolipids as non viral vectors for DNA transfection in hematopoietic cell lines and CD34+ cells. *Blood Cells Molec. & Diseases* 23:69-87.
2. Guillaume-Gable *et al.*, 1998. Cationic phosphonolipids as nonviral gene transfer agents in the lung of mice. *Hum Gene Ther* 9:2309-19.
3. Xie M. *et al.*, 2016. PKM2-dependent glycolysis promotes NLRP3 and AIM2 inflammasome activation. *Nat Commun.* 7:13280.
4. Järver P. *et al.*, 2018. Single-Stranded Nucleic Acids Regulate TLR3/4/7 Activation through Interference with Clathrin-Mediated Endocytosis. *Sci Rep.* 8:15841.
5. Katashiba Y. *et al.*, 2011. Interferon-α and interleukin-12 are induced, respectively, by double-stranded DNA and single-stranded RNA in human myeloid dendritic cells. *Immunology.* 132(2):165-73.

TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873

InvivoGen USA (International): +1 (858) 457-5873

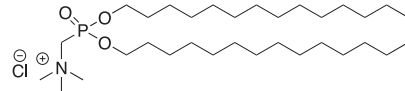
InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Asia: +852 3622-3480

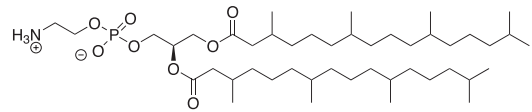
E-mail: info@invivogen.com

STRUCTURE

LyoVec™ is a combination of the phosphonolipid DTCPTA and the neutral lipid DiPPE (structures below).



Di-tetradecylphosphoryl-N,N,N-trimethylmethanaminium chloride (DTCPTA)



1,2-Diphytanoyl-sn-Glycero-3-Phosphoethanolamine (DiPPE)

APPLICATIONS

LyoVec™ can be used for adherent cells as well as cells grown in suspension. It has been validated for use as a:

- Cationic lipid-based transfection reagent
- Nucleic acid complexing agent

METHODS

Summary for preparing LyoVec™-nucleic acid complexes



01

Reconstitute lyophilized LyoVec™ with sterile H₂O and leave for 30 min at 15-25°C

02

Prepare the complexes at 10 µg/ml

- Add 1 µl of nucleic acid (1 mg/ml) to a 1.5 ml tube
- Add 100 µl of LyoVec™
- Mix gently by tapping the tube
- Incubate for 15 min -1 hour at 15-25°C
- Use immediately



LyoVec™ reconstitution

- Resuspend with deionized sterile water
 - Add 2 ml per small vial (cat. codes: **lyec-1** and **lyec-2**).
 - Add 8 ml per large vial (cat. code: **lyec-3**).
- Homogenize by vortexing gently for 30 seconds.
- Leave vial at room temperature (15-25°C) for at least 30 minutes prior to initial use. Use immediately or store at 4°C.

The protocol for the preparation of LyoVec™-nucleic acid complexes and for the transfection of mammalian cells is given on the next page.

Protocol for the preparation of LyoVec™-nucleic acid complexes

1. If stored at 4°C, bring LyoVec™ to room temperature and gently vortex to homogenize before use.
 2. In a sterile 1.5 ml microfuge tube at room temperature, add 1 µl of nucleic acid (1 mg/ml). Add 100 µl of LyoVec™. Mix gently by tapping the tube.
 3. Incubate at room temperature (15-25°C) for 15 minutes to 1 hour to allow for the formation of the complexes.
 4. Use the LyoVec™-nucleic acid complexes immediately.
- Note:* Do not keep LyoVec™-nucleic acid complexes for more than 2 hours.

Transfection of cells

The table below shows indicative values for the volume of medium, the volume of LyoVec™/DNA complex and the number of cells for common transfection experiments.

Culture format	96-well plate	24-well plate	12-well plate	6-well plate	100 mm dish
Medium per well/dish	200 µl	500 µl	1 ml	2 ml	8 ml
LyoVec™/DNA complex	10 µl	25 µl	50 µl	100 µl	400 µl
Cells per well/dish	1-5 x 10 ⁴	1-2 x 10 ⁵	2-4 x 10 ⁵	5-10x10 ⁵	3-5 x 10 ⁶

1. Seed cells in culture medium containing serum in the dish or well as specified in the table above. Cells should be 50-80% confluent for best results. The transfection can be performed immediately after cell seeding.

Note: Cells can be added to dish or well before or after LyoVec™/DNA complex without affecting transfection efficiency.

2. Add LyoVec™-DNA complexes drop by drop to the medium at a 1/20 volume ratio (see table above). Swirl to distribute the complexes uniformly.
3. Incubate under appropriate conditions for 24-72 hours to allow for gene expression then assay or apply antibiotic for selection.

For optimal transfection:

The ratio volume of cells/volume of complex is an important parameter for achieving optimal transfection, and may need to be optimized for your cells. In optimization experiments, test various cell/complex ratios by reducing or increasing the volume of complexes. Also note that LyoVec™ complexes work best at ratios for which a slight toxicity is observed.

RELATED PRODUCTS

Product	Description	Cat. Code
Cytosolic DNA sensors (CDSs) ligands		
G3-YSD	Y-form DNA	ttrl-ydna
HSV-60 Naked	Viral DNA motif	ttrl-hsv60n
VACV-70 Naked	Viral DNA motif	ttrl-vav70n
Endosomal TLR ligands		
ssPoly(U) Naked	TLR7/TLR8 ligand	ttrl-sspu
RIG-I ligands		
3p-hpRNA	5' triphosphate hairpin RNA	ttrl-hprna
5'-ppp-dsRNA	5' triphosphate dsRNA	ttrl-3prna
Poly(dA:dT)	Synthetic analog of B-DNA	ttrl-patn
Poly(I:C) HMW	Synthetic dsRNA polymer	ttrl-pic
Expression plasmids		
pUNO1-mcs	Cloning plasmid	puno1-mcs
pMONO-zeo-mcs	For a single gene	pmonoz-mcs
pDUO-mcs	For two genes	pduo-mcs
Expression plasmids - to generate tagged proteins		
pSELECT-NGFP-zeo	GFP-N-terminal tag	psetz-ngfp
pSELECT-NHA-zeo	HA-N-terminal tag	psetz-nha
pSELECT-NHis-zeo	HIS-N-terminal tag	psetz-nhis
pSELECT-NLucia-zeo	Lucia-N-terminal tag	psetz-nlucia
Expression plasmids - for antibody production		
pFUSE-CHlg-hG1	Heavy chain, human IgG1	pfuse-hchg1
pFUSE2-CLlg-hK	Light chain, human Kappa	pfuse2-hclk
pTRIOZ-hlgG1	Heavy & light chains	p trio z-higg1

For a complete list of our cloning and expression vectors, visit <https://www.invivogen.com/vectors>

TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873
InvivoGen USA (International): +1 (858) 457-5873
InvivoGen Europe: +33 (0) 5-62-71-69-39
InvivoGen Asia: +852 3622-3480
E-mail: info@invivogen.com