



# EndoFectin™ Transfection Reagents

EndoFectin™ Max is a lipid-based transfection reagent that has been optimized for highly efficient transfection of a large collection of commonly used cell lines as well as many difficult-to-transfect cell lines.

## Advantages

- ◆ Superior transfection efficiency
- ◆ Easy to use
- ◆ Low cytotoxicity



## Versatile & Efficient For Robust Transfection

- √ Effectively deliver plasmid, siRNA (gene knockdown) and Cas9/sgRNA ribonucleoprotein particles (RNPs) to many cell types.
- √ Better transfection efficiency than LP2000 and LP3000 for a broad range of cell lines.

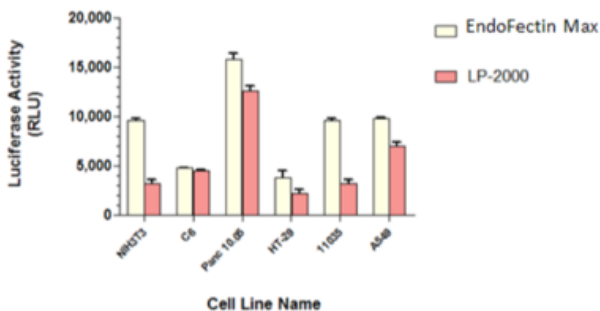
## Easy To Use

Does not require removal of serum or culture medium.

## Low Cytotoxicity

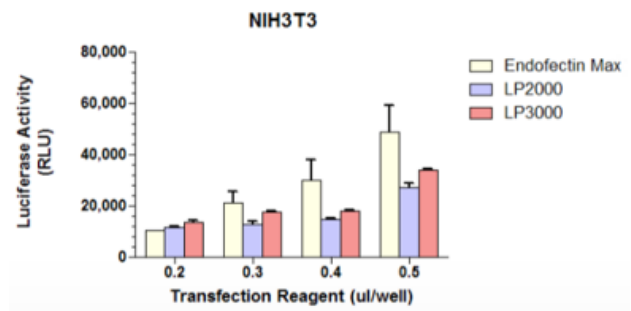
Retain more live cells after transfection.

**Comparison of EndoFectin™ Max and Lipofectamine® 2000 in plasmid transfection**



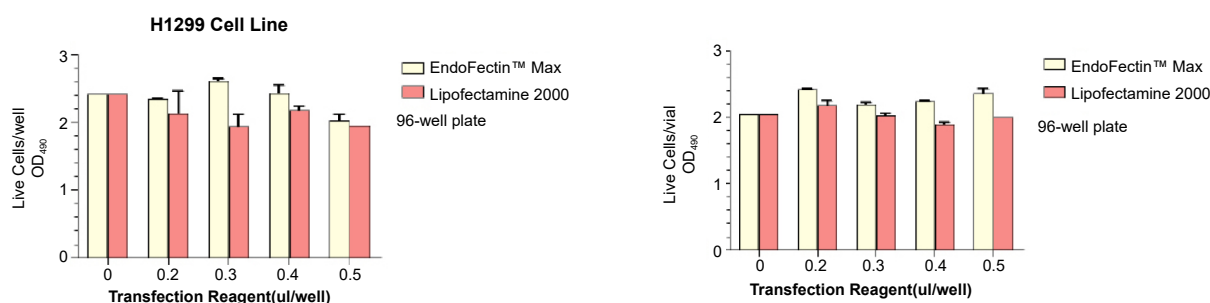
**Figure 1.** Transfection of 7 commonly-used cell lines using either EndoFectin™ Max or Lipofectamine® 2000 (LP-2000). This image shows cell lines with high transfection efficiency and were transfected with a firefly luciferase-expressing plasmid. Transfection efficiencies were assessed by luciferase activity (RLU). Lipofectamine® is a trademark of Invitrogen.

**Comparison of EndoFectin™ Max and Lipofectamine® 3000 in plasmid transfection**



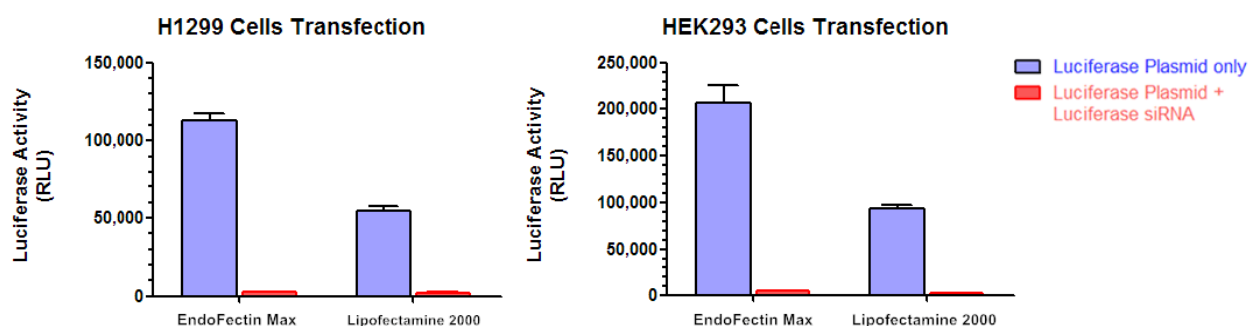
**Figure 2.** Transfection of NIH3T3 cells using either EndoFectin™ Max, Lipofectamine® 2000 (LP2000) or Lipofectamine® 3000 (LP3000). NIH3T3 was transfected with a firefly luciferase-expressing plasmid. Transfection efficiencies were assessed by luciferase activity (RLU).

## Comparison of EndoFectin™ Max and Lipofectamine® 2000 for cytotoxicity



**Figure 3.** Comparison of cytotoxicity between EndoFectin™ Max and Lipofectamine® 2000. Lipofectamine® is a trademark of Invitrogen.

## Comparison of EndoFectin™ Max and Lipofectamine® 2000 in siRNA transfection



**Figure 4.** Transfection of two commonly-used cell lines (H1299 & HEK293) using EndoFectin™ Max or Lipofectamine® 2000. Cell lines were transfected with: a. A firefly luciferase-expressing plasmid (blue bars), b. Firefly luciferase-expressing plasmid and Luciferase siRNA co-transfected (red bars). Transfection efficiencies were assessed by luciferase activity (RLU). Lipofectamine® is a trademark of Invitrogen.

Endofectin Products	Application	Catalog number	Price
EndoFectin™ Max	High-efficiency transfection over a broad range of cell lines	EF013 (1 mL)	\$295
		EF014 (3 mL)	\$589
EndoFectin™ Expi293	Optimized for transfection of the suspension cell line Expi293F	EF007 (1 mL)	\$320
		EF009 (1L culture)	\$485
		EF010 (10L culture)	\$4155
EndoFectin™ HepG2	Optimized for transfection of the hepatocellular carcinoma cell line HepG2	EF005 (1 mL)	\$355
		EF006 (3 mL)	\$705
EndoFectin™ Lenti	Transfection of cells for high-titer lentiviral packaging	EF001 (1 mL)	\$245
		EF002 (3 mL)	\$489

### Contact us:

#### GeneCopoeia, Inc.

9620 Medical Center Dr. Suite 101,  
Rockville, MD 20850 USA



Email: [inquiry@genecopoeia.com](mailto:inquiry@genecopoeia.com)  
Tel: +1 (301) 762-0888  
Fax: +1 (301) 762-3888  
Website: [www.genecopoeia.com](http://www.genecopoeia.com)