



EndoFectin™ Max Transfection Reagents

Superior Transfection Efficiency

Superior transfection efficiency in both DNA and siRNA transfection.
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.

Superior Transfection efficiency in many difficult-to-transfect cell lines

Greater transfection performance in 6 hard-to-transfect cell lines.

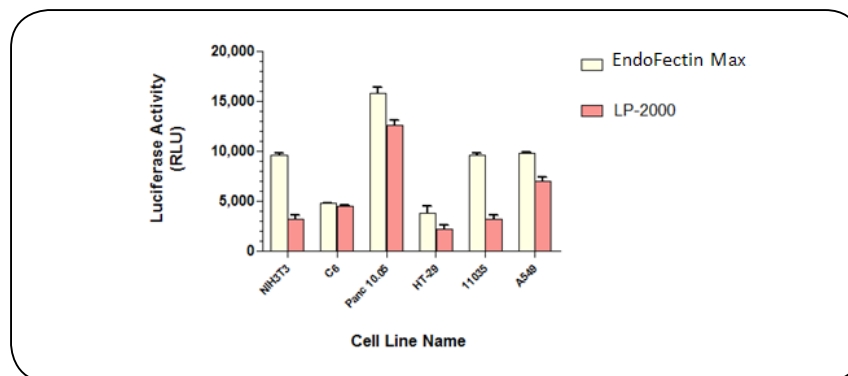


Figure 1. Transfection of 6 commonly-used cell lines using either EndoFectin™ Max or Lipofectamine®2000 (LP-2000). Cells following transfection with a firefly luciferase-expressing plasmid were analyzed for luciferase activity (RLU). Lipofectamine® is a trademark of Invitrogen.

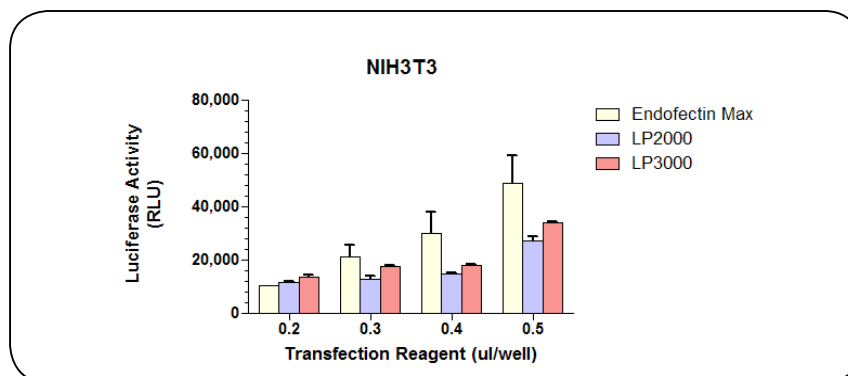


Figure 2. Transfection of NIH3T3 cells using either EndoFectin™ Max, Lipofectamine® 2000 (LP2000) or Lipofectamine® 3000 (LP3000). NIH3T3 was transfected with a firefly luciferase-expressing plasmid. Transfection efficiencies were assessed by luciferase activity (RLU).

Superior efficiency in siRNA transfection

Higher transfection efficiency compared with Lipofectamine® 2000 in siRNA transfection.

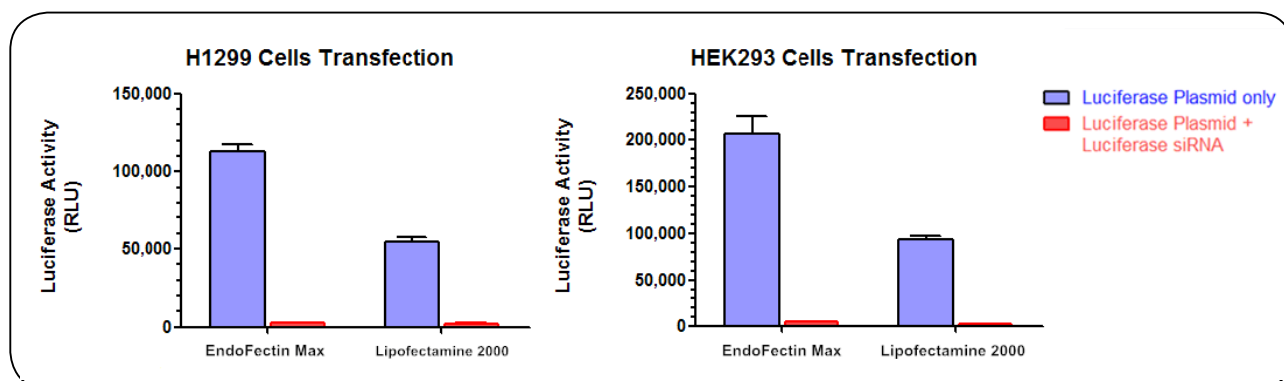


Figure 3. Transfection of two commonly-used cell lines (H1299 & HEK293) using EndoFectin™ Max or Lipofectamine® 2000. Cell lines were transfected with: a. A firefly luciferase-expressing plasmid (blue bars), b. Firefly luciferase-expressing plasmid and Luciferase siRNA co-transfected (red bars). Transfection efficiencies were assessed by luciferase activity (RLU). Lipofectamine® is a trademark of Invitrogen.

Low cytotoxicity

Higher cell viability after transfection.

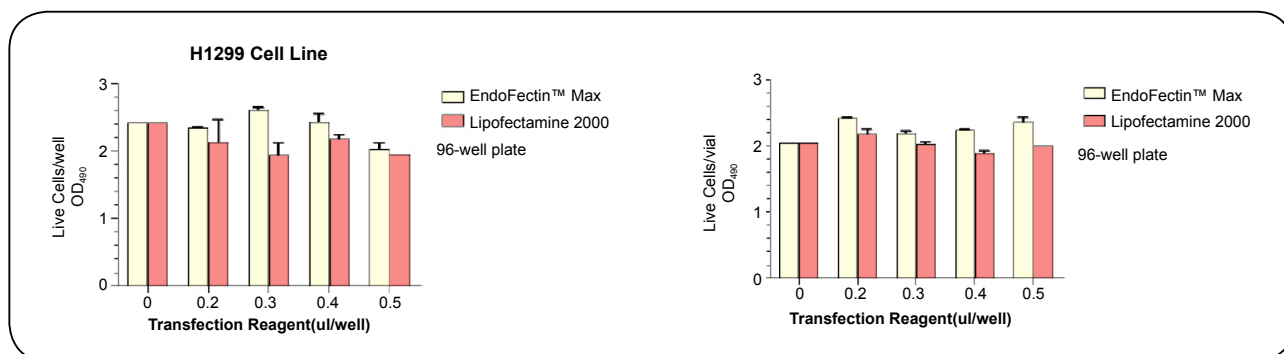


Figure 4. Comparison of cytotoxicity between EndoFectin™ Max and Lipofectamine® 2000. Lipofectamine® is a trademark of Invitrogen.

Endofectin Products	Application	Catalog number	Price
EndoFectin™ Max	High-efficiency transfection over a broad range of cell lines	EF003 (1 mL)	\$295
		EF004 (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

To Order

Visit our website.

<http://www.genecopoeia.com/product/endofectin/>

For more information please contact us at

inquiry@genecopoeia.com

or 1-866-360-9531

EF-031318