





## Lipid Transfection Reagents Selection Guide

Bio-Rad offers a family of lipid transfection reagents that effectively deliver nucleic acids to a wide range of cultured mammalian cell lines, including ones that are difficult to transfect. Each lipid is optimal for certain applications. The siLentFect lipid reagent is an RNAi-specific lipid that is ideal for siRNA delivery. It can also be used to perform co-transfection with a small reporter-plasmid along with the siRNA molecule. The TransFectin lipid reagent is a general-purpose lipid that is ideal for plasmid delivery in a wide range of cell lines. The cell line-specific lipid reagents, COSFectin and HEKfectin, are best suited for experiments using the specified cell line and are particularly valuable for protein over-expression and viral production.

	siLentFect Lipid Reagent	TransFectin Lipid Reagent	HEKfectin Cell Line-Specific Lipid Reagent	COS Fectin Cell Line-Specific Lipid Reagent
				
Purpose	RNAi-specific lipid	General-purpose lipid	HEK cell line-specific lipid	COS cell line-specific lipid
Applications	RNAi (siRNA delivery)*	Gene expression RNAi (siRNA/shRNA delivery)	Protein over-expression Viral production**	Protein over-expression Viral production**
Cell lines***	Variety of adherent and suspension cell lines, including A549, BHK, HeLa, Jurkat, K562, primary fibroblasts, and Vero	Variety of adherent and suspension cell lines, including A549, HeLa, HUVEC, Jurkat, MCF-7, and primary fibroblasts	HEK 293 cells and subclones (H, T, F, and Phoenix)	COS cells and subclones (1, 3, and 7)
Shelf life	6 months	6 months	6 months	6 months

\* siLentFect lipid reagent can successfully co-transfect an siRNA molecule with a plasmid DNA.

\*\* Not for gene expression.

\*\*\* For a more complete list of cell lines, go to [www.bio-rad.com/transfection/](http://www.bio-rad.com/transfection/).



**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

---

**Web site** [www.bio-rad.com](http://www.bio-rad.com) **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 11 5044 5699  
**Canada** 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00  
**France** 01 47 95 69 65 **Germany** 089 31 884 0 **Greece** 30 210 9532 220 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300  
**Israel** 03 963 6050 **Italy** 39 02 216091 **Japan** 03 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666  
**New Zealand** 64 9 415 2280 **Norway** 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04  
**Singapore** 65 6415 3188 **South Africa** 27 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 061 717 95 55  
**Taiwan** 886 2 2578 7189 **Thailand** 800 88 22 88 **United Kingdom** 020 8328 2000

---