



# Powerful Delivery

TransFectin™ Lipid Reagent



High efficiency, low cytotoxicity, and a simple protocol put Bio-Rad's TransFectin lipid reagent in a league of its own.





# TransFectin lipid reagent for powerful, efficient transfection

For efficient and reproducible transfection, Bio-Rad offers TransFectin, a powerful cationic lipid reagent effective in difficult-to-transfect cell lines as well as in primary cells. TransFectin reagent's advanced formulation delivers the highest efficiencies and transgene activity in a broad range of cell lines while minimizing damaging cytotoxic effects observed with other high-efficiency transfection reagents.

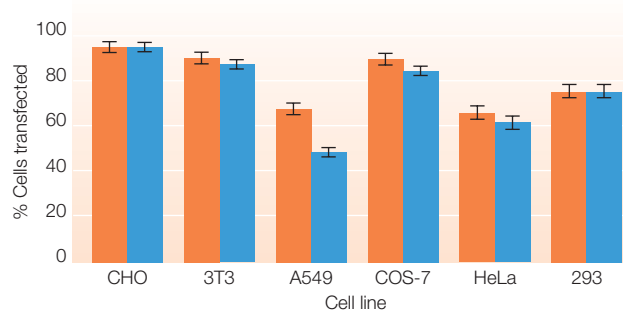
## Features

- High efficiency and activity for both easy- and difficult-to-transfect cells
- Lower cytotoxicity — maintains cell viability
- Excellent performance at culture densities between 40 and 90%
- Simple 3-step protocol — no media changes required for most cell types

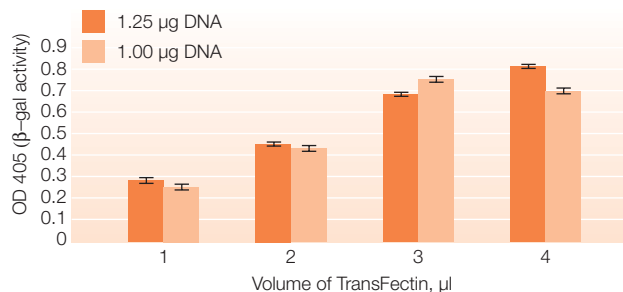
TransFectin reagent allows high transfection efficiencies with lower amounts of lipid than most other lipid reagents. This means more experiments per vial of reagent and lower costs per transfection.

## Ordering Information

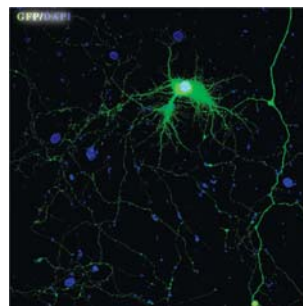
Catalog #	Description
170-3350	TransFectin Lipid Reagent, 0.5 ml
170-3351	TransFectin Lipid Reagent, 1.0 ml
170-3352	TransFectin Lipid Reagent, 5 x 1.0 ml



**Comparison of transfection efficiency using TransFectin (■) or another popular lipid transfection reagent (■).** Cell lines were transfected with pCMV.SPORT-β-gal and stained with X-gal 24 hr after transfection. Efficiency is reported as the percentage of cells in the population expressing β-galactosidase.



**Effective delivery to suspension cells using TransFectin.** K562 cells were plated in 24-well plates at a density of  $4 \times 10^5$  cells/well. TransFectin:plasmid DNA complexes were formed by combining 1.25 μg (■) or 1.00 μg (■) pCMV.SPORT-β-gal with TransFectin. After 20 min, transfection complexes were added to the cells in the presence of serum. At ~24 hr posttransfection, cells were pelleted and lysed, then assayed for β-gal expression.



## Efficient transfection of primary neuronal cells with TransFectin.

Primary rat hippocampal neurons were transfected on coverslips in 24-well plates using 1 μg of a GFP expression plasmid and 2 μl TransFectin. At 3 hr after complex addition, the medium was replaced with 50:50 conditioned: fresh medium. Neurons were fixed and analyzed 48 hr after transfection. Data courtesy of Dr Jamil Kanaani, Diabetes Center, University of California, San Francisco.

## Examples of Cell Lines Used

104C1
143B
A-498
A-549
BHK
CHO
CHO-DG44
COS-7
CV-1
DT40
F9
HCT 116
HEK 293
HeLa
Hs68—primary human fibroblasts
HT-22
HT-1080
Jurkat
K-562
K-B
LLC-MK2
LNCaP
MCF-7
MDCK
MIN6
Neuro-2a
NIH 3T3
PC-3
PC-12
Breast—primary human fibroblasts
Primary rat myoblasts
Primary rat hippocampal neurons
SH-SY5Y
SK-HEP-1
U-2 OS
U-87 MG
Vero



**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Web site [www.bio-rad.com](http://www.bio-rad.com) USA 800 4BIORAD Australia 02 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 21 3237 9400  
Canada 905 712 2771 China 86 21 6426 0808 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65  
Germany 089 318 84 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 455 8800 India 91 124 4029300/5013478 Israel 03 963 6050  
Italy 39 02 216091 Japan 03 5811 6270 Korea 82 2 3473 4460 Mexico 55 5200 05 20 The Netherlands 0318 540666 New Zealand 64 9415 2280  
Norway 23 38 41 30 Poland 48 22 331 99 99 Portugal 351 21 472 7700 Russia 7 095 721 14 04 Singapore 65 6415 3188 South Africa 27 0861 246 723  
Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 061 717 95 55 Taiwan 886 2 2578 7189/2578 7241 United Kingdom 020 8328 2000