



TransFectin™ Lipid Reagent Protocols

Cell type:	A549
Species:	Human
Tissue:	Lung; carcinoma
Primary or established:	Established
Adherent or suspension:	Adherent

Nucleic Acid Characteristics

Nucleic acid type:	Plasmid	Nucleic acid isolation method:	Aurum plasmid kit
Vector:	pCMV Sport β -Gal		

Growth Conditions

Vessel type and size:	24-well plate	Growth medium:	D-MEM, 0.5 ml
Number of cells seeded:	$2.5\text{--}5.0 \times 10^4$	Serum concentration:	10%
Growth period prior to transfection:	24 hr		

Transfection Conditions

Cell density at transfection:	50–70%	Volume of TransFectin used (per well):	0.5–1.0 μ l
Amount of DNA used (per well or vessel):	0.5–1.0 μ g	Length of complex incubation with cells:	24 hr

Analysis

Time between transfection and analysis:	24 hr	Transfection efficiency:	60–70%
Reporter assay used:	β -Gal, ONPG		

Protocol Notes

Protocol number:	15	Author's name:	Sean Baker
		Institution:	Bio-Rad Laboratories