



TransFectin™ Lipid Reagent Protocols

Cell type:	PC12
Species:	Rat
Tissue:	Pheochromocytoma
Primary or established:	Established
Adherent or suspension:	Adherent

Nucleic Acid Characteristics

Nucleic acid type:	Plasmid DNA	Nucleic acid isolation method:	Not specified
Vector:	pEGFP-C1		

Growth Conditions

Vessel type and size:	12-well plate	Growth medium:	D-MEM
Number of cells seeded:	1.5 x 10 ⁵	Serum concentration:	15%
Growth period prior to transfection:	24 hr		

Transfection Conditions

Cell density at transfection:	70–90%	Volume of TransFectin used (per well):	2 µl
Amount of DNA used (per well or vessel):	1 µg	Length of complex incubation with cells:	24 hr

Analysis

Time between transfection and analysis:	48 hr	Transfection efficiency:	40–50%
Reporter assay used:	GFP fluorescence		

Protocol Notes

Protocol number:	5	Author's name:	Anonymous
		Institution:	