



TransFectin™ Lipid Reagent Protocols

Cell type:	143B
Species:	Human
Tissue:	Bone marrow, osteosarcoma
Primary or established:	Established
Adherent or suspension:	Adherent

Nucleic Acid Characteristics

Nucleic acid type:	Plasmid	Nucleic acid isolation method:	Promega Wizard kit
Vector:	pEGFP, pDsRED		

Growth Conditions

Vessel type and size:	6-well plate	Growth medium:	D-MEM
Number of cells seeded:	2.0–5.0 x 10 ⁵	Serum concentration:	10%
Growth period prior to transfection:	24–48 hr		

Transfection Conditions

Cell density at transfection:	50–60%	Volume of TransFectin used (per well):	2–5 µl
Amount of DNA used (per well or vessel):	1 µg	Length of complex incubation with cells:	3–24 hr

Analysis

Time between transfection and analysis:	24–48 hr	Transfection efficiency:	60–90%
Reporter assay used:	GFP or DsRED fluorescence		

Protocol Notes Removal of complexes after 3 hr reduces cytotoxicity.

Protocol number:	9	Author's name:	Anonymous
		Institution:	FinMIT Center of Excellence