



## TransFectin™ Lipid Reagent Protocols

<b>Cell type:</b>	HEK 293T
<b>Species:</b>	Human
<b>Tissue:</b>	Kidney
<b>Primary or established:</b>	Established
<b>Adherent or suspension:</b>	Adherent

---

### Nucleic Acid Characteristics

<b>Nucleic acid type:</b>	Plasmid	<b>Nucleic acid isolation method:</b>	Not specified
<b>Vector:</b>	pEGFP-N1		

---

### Growth Conditions

<b>Vessel type and size:</b>	24-well plate	<b>Growth medium:</b>	D-MEM
<b>Number of cells seeded:</b>	9.0 x 10 <sup>4</sup> cells/well	<b>Serum concentration:</b>	10%
<b>Growth period prior to transfection:</b>	24 hr		

---

### Transfection Conditions

<b>Cell density at transfection:</b>	80%	<b>Volume of TransFectin used (per well):</b>	0.5–1.25 µl
<b>Amount of DNA used (per well or vessel):</b>	0.25–0.5 µg	<b>Length of complex incubation with cells:</b>	12 hr

---

### Analysis

<b>Time between transfection and analysis:</b>	54 hr	<b>Transfection efficiency:</b>	80–90%
<b>Reporter assay used:</b>	GFP		

---

### Protocol Notes

---

<b>Protocol number:</b>	17	<b>Author's name:</b>	Anonymous
		<b>Institution:</b>	Ein-Karem Medical School