

Polycystin1 the Gene Product of PKD1 Induces Resistance to Apoptosis and Spontaneous...

Polycystin1 the Gene Product of PKD1 Induces Resistance to Apoptosis and Spontaneous Tubulogenesis in MDCK Cells

ABSTRACT

The authors use the DeadEnd™ Fluorometric TUNEL System (Cat.# G3250) to examine the effect of expression of PKD1 in MDCK cells on the induction of apoptosis.

Cell Notes 2, 10.

Publication Date: 2000



Download Article

(126 KB)