

# An Alternative NonApoptotic Form of Programmed Cell Death

---

## ABSTRACT

*The authors use DeadEnd™ Fluorometric TUNEL System (Cat.# G3250) to investigate events associated with IGFIR-IC-induced cell death. IGFIR-IC is human insulin-like growth factor I receptor beta subunit intracellular domain.*

*Cell Notes* **2**, 9.

**Sperandio, S. et al. (2000), 14376–14381 Review by Rebecca Sanders\***

*Proc. Natl. Acad. Sci. USA* **20**, Agnes Scott College, Decatur GA, USA.\*

Publication Date: 2000



## Download Article

(89 KB)

---