



**Dr. Stewart Abbot**  
**Senior Director Stem Cells, Celgene Cellular Therapeutics (USA)**

### **Biography**

*Dr. Stewart Abbot* is Senior Director of Stem Cell Research (Celgene Cellular Therapeutics). He holds a B.Sc. in Biological Sciences (Edinburgh), M.Sc. in Biomedical Engineering (Glasgow) and Ph.D. in Pathology (London). His academic career focused on vascular biology, pharmacology and toxicology. He joined Amersham Biosciences (GE Healthcare) in 2000 to develop cell-based assay systems for pharmaceutical screening. Following acquisition of Amersham he transferred to General Electric's Global Research Center (Albany, NY) to manage the Molecular and Cellular Biology research laboratory. In this laboratory Dr. Abbot developed GE's expertise in umbilical cord blood and human embryonic stem cell biology. In 2007 he joined Celgene Cellular Therapeutics and now leads the group's efforts to isolate, characterize and develop novel therapeutics based on human placenta-derived stem cells. This work is expected to result in clinical trials early in 2009. Throughout his career he has worked across the traditional boundaries of life and physical sciences and has managed a number of successful basic science and cell-based product development programs.