



**Dr. Davinder Gill**  
**Assistant Vice President, Biologic Therapeutics, Wyeth (USA)**

### **Biography**

Davinder Gill is Assistant Vice President, Biologic Therapeutics at Wyeth Research. His group is responsible for developing protein therapeutics within Wyeth's Discovery pipeline. Since joining in 2003, Davinder's group has advanced 15 proteins into Development, 7 of which are in various stages of human clinical trials. Prior to joining Wyeth, Davinder was Associate Director, Biotherapeutic Discovery at Millennium Pharmaceuticals in Cambridge, Massachusetts. Davinder carried out a postdoctoral fellowship at Harvard Medical School. Davinder is a Chemical Engineer by training. He obtained a PhD from the University of Houston, Texas and a Bachelors degree from the Indian Institute of Technology (IIT).