

**Characterisation of the 5-HT<sub>6</sub> receptor using CatchPoint  
cAMP assay: comparison with FP and FlashPlate.**

Martyn Wood, Lindsay Mulligan and Nisha Patel.

GlaxoSmithKline,

New Frontiers Science Park, Third Avenue, Harlow, Essex, CM19  
5AW, U.K.

Traditional methods for the study of GS and GI coupled receptors have followed changes in the level of cAMP using radioactive methods, either pre-labeling using <sup>3</sup>H-adenine or using antibody-competition with labeled conjugate e.g. <sup>125</sup>I-cAMP. Recently, new methodologies have been developed which allow the measurement of cAMP levels using non-radiometric determination. These methods include fluorescence, fluorescence polarisation (FP) and luminescence proximity. We have therefore evaluated these methodologies in order to develop a non-radioactive medium throughput assay in intact cells. We have used NEN FlashPlate and NEN FP techniques and compared them to Molecular Devices CatchPoint ELISA assay, using the human 5-HT<sub>6</sub> receptor expressed in HeLa cells. We have found the CatchPoint assay to be a sensitive and robust assay which offers many advantages.