

## Development of a Rapid Capture ELISA Using PCR Products and the PinPoint™ System

## **ABSTRACT**

A simple procedure for the development of an ELISA is presented. The procedure uses the PinPoint™ Xa-1 T-Vector System to express a protein antigen encoded by a PCR product. Because the PinPoint™ System adds a biotin tag to the protein, a single-step purification by affinity for streptavidin allows direct use in ELISA. We report the use of this procedure to express a recombinant protein from the nucleocapsid domain of the feline foamy virus (FeFV) gag gene, fused with a biotin tag. This fusion protein was applied directly to streptavidin-coated ELISA wells. Antibody to FeFV was detected in this ELISA with a 100% correlation to other detection methods, including immunoblot, serum neutralization and virus isolation.

## Promega Notes 70, 14.

Ingrid G. Winkler1, Jean-Pierre Levesque2 and Robert L.P. Flower3

1School of Pharmacy and Medical Science, University of South Australia; 2Haematology Division, Hanson Centre for Cancer Research, Australia; 3Pacific Laboratory Medicine Services, Royal North Shore Hospital, Australia
Publication Date: 1999



## **Download Article**

(279 KB)