

SignaTECT™ DNA Dependent Protein Kinase Assay System

ABSTRACT

The SignaTECT™ DNA-Dependent Protein Kinase Assay System developed at Promega accurately measures the protein kinase activity of DNA-dependent protein kinase (DNA-PK), both in purified samples and nuclear extracts. The assay can be performed quickly and gives very low background counts indicating minimal interference from endogenously phosphorylated proteins present in nuclear extracts. This new assay system provides the researcher a powerful tool with which to measure DNA-PK activity.

Promega Notes 63, 7.

Said Goueli and Kevin Hsiao

Promega Corporation

Publication Date: 1997



Download Article

(179 KB)
