The Role of Cell-Free Rabbit Reticulocyte Expression Systems in Functional Genomics

Article

Comments & Ratings

ABSTRACT

In the broad area of functional genomics, that is the global characterization of proteins and their function, cell-free rabbit reticulocyte lysate (*RRL*) has been used extensively to elucidate the mechanisms of mammalian translation, cotranslational modifications, post-translational modifications and translocation of proteins. More recently, *RRL* has been used as the workhorse for manufacturing the proteins engaged in interaction, selection and protein evolution studies from DNA or mRNA libraries either in microarray, display or in vitro expression cloning (*IVEC*) technologies. This chapter highlights recent functional proteomics applications that use cell-free mammalian *RRL*.

Michele Arduengo, Elaine Schenborn and Robin Hurst Promega Corporation



Download Article

(1 MB)



It appears that you have Javascript disabled. Our website requires Javascript to function correctly. For the best browsing experience, please enable Javascript.