

Bioluminescence & Chemiluminescence

Molecular reporting from photons

Ed.: J.W.Hastings L.J.Kricka P.E.Stanley, Chichester: Wiley 1996: 509-512

Chemiluminescence in situ hybridization for the detection of viral genomes

Pasini P, M Musiani, A Roda, M Zerbini, G Gentilomi, M Baraldini, G Gallinella and S Venturoli
University of Bologna

Keywords: LB 980 Luminograph, chemiluminescence in situ hybridization, Parvovirus, bone marrow cells