

Detection of Dipeptidyl Peptidase Activity with DPPIV-Glo™ Assay

ABSTRACT

Ranbaxy Laboratories Ltd., is a research-based, international pharmaceutical company located in India. In addition to being a dominant player in the area of developing generic medicines, Ranbaxy has committed to investing in drug discovery. Keeping in mind the highly complex, cost-intensive and failure-prone nature of drug discovery, Ranbaxy has evolved an alliance-centric business model to bring sustainability to its drug discovery effort. Ranbaxy has entered into alliance with partners at different stages of drug discovery and development. A partnership with Promega is an effort to access robust assays that can assist Ranbaxy in hit identification and lead optimization during early drug discovery stages. Here we describe our characterization of the DPPIV-GLOTM Protease Assay.

Cell Notes 23, 6-9.

Joseph Alex Davis, Vamshi Krishna Tulasi Abhijit Ray, and Pradip Bhatnagar Ranbaxy Laboratories, Ltd. India
Publication Date: 2009



Download Article

(301 KB)

HOW TO CITE THIS ARTICLE

Scientific Style and Format, 7th edition, 2006

Davis, J. A., Ray, V. K. T. A and Bhatnagar, P. Detection of dipeptidyl peptidase activity with DPPIV-Glo™ Assay. [Internet] 2009. [cited: year, month, date]. Available from: http://www.promega.com/resources/pubhub/cellnotes/detection-of-dipeptidyl-peptidase-activity-with-dppiv-glo-assay/

American Medical Association, Manual of Style, 10th edition, 2007

Davis, J. A., Ray, V. K. T. A and Bhatnagar, P. Detection of dipeptidyl peptidase activity with DPPIV-Glo™ Assay. Promega Corporation Web site. http://www.promega.com/resources/pubhub/cellnotes/detection-of-dipeptidyl-peptidase-activity-with-dppiv-glo-assay/ Updated 2009. Accessed Month Day, Year.