

DDP-IX Protease Assay Service

Dipeptidyl peptidase IX [Protease Assay](#) for compound screening and profiling via fluorescence-based quantification of enzyme activity.

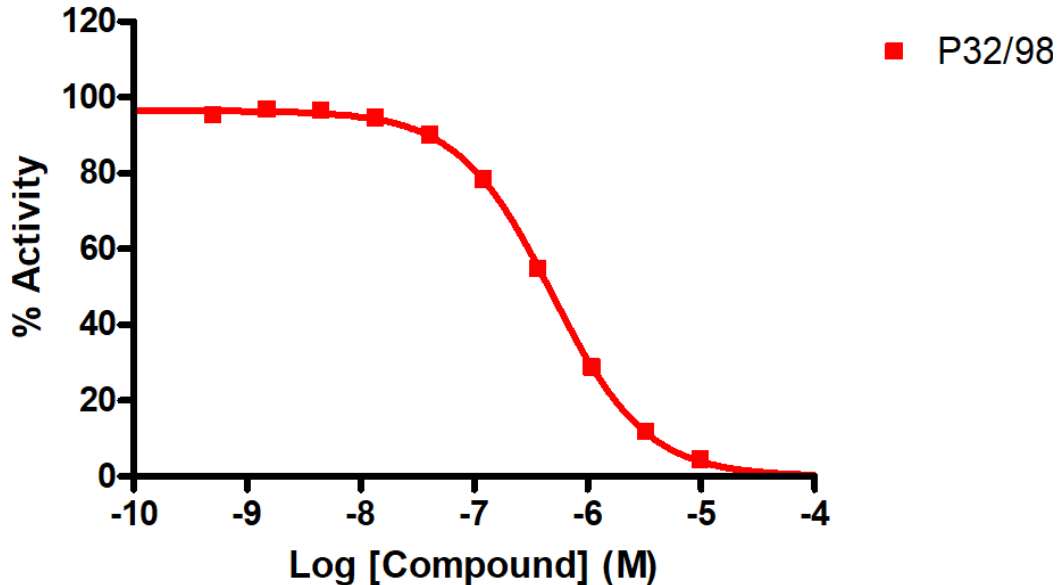
Target Overview

CAT#	DPP-IX
Enzyme	Human Dipeptidyl peptidase IX
Synonym	Dipeptidyl peptidase IX, DPP9, CD26, DPLP9, DPRP2
Substrates	H-GP-AMC, 10 μ M

Assay Properties

Readout	Enzymatic activity
Reference Compound IC50s	Compounds IC50 (nM) P32-98 700
Scientific Information	Q86T12 (https://www.uniprot.org/uniprotkb/Q86T12/entry)
Screening Location	Malvern, PA, USA
Further information	More information can be found on our website Protease Assay Services .

Reference compound IC50 for DPP-IX



Let's Discover Together.

© Jun 2023 **Reaction Biology**. All rights reserved.

https://www.reactionbiology.com/datasheet/dpp-ix_protea_malvern

