

DPP-VIII Protease Assay Service

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

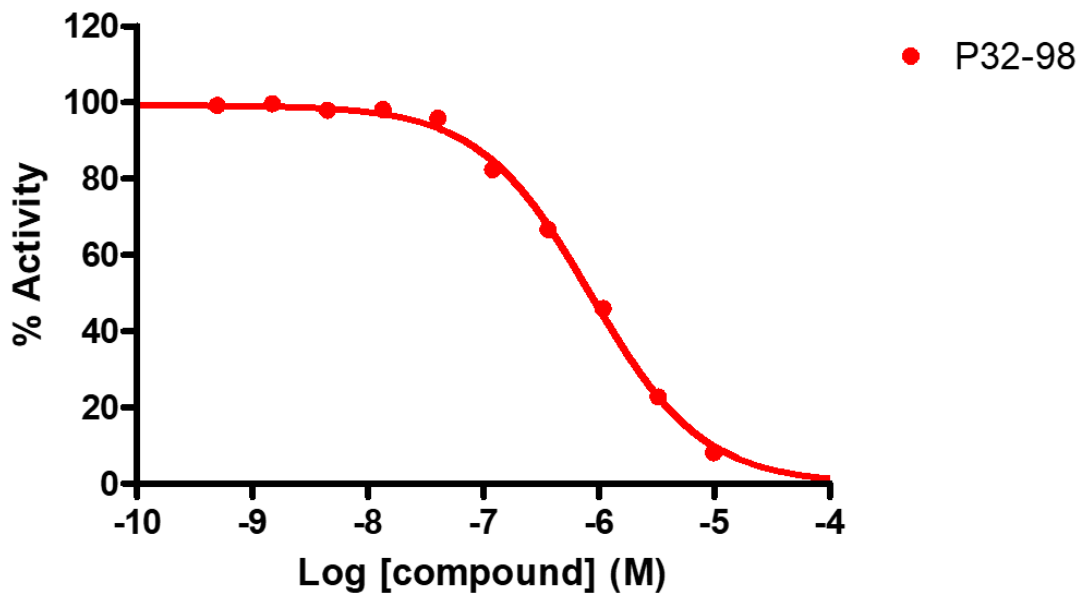
Target Overview

CAT#	DPP-VIII
Enzyme	Human Dipeptidyl peptidase VIII
Synonym	Dipeptidyl peptidase 8, Dipeptidyl peptidase IV-related protein 1, MSTP097, MSTP135, MSTP141
Substrates	H-GP-AMC, 10 μ M

Assay Properties

Readout	Enzymatic activity
Reference Compound IC50s	Compounds IC50 (nM) P32-98 967
Scientific Information	Q6V1X1 (https://www.uniprot.org/uniprotkb/Q6V1X1/entry)
Screening Location	Malvern, PA, USA
Further information	More information can be found on our website Protease Assay Services .

Reference compound IC50 for DPP-VIII



Let's Discover Together.

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

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

