

# DPP-IV Protease Assay Service

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

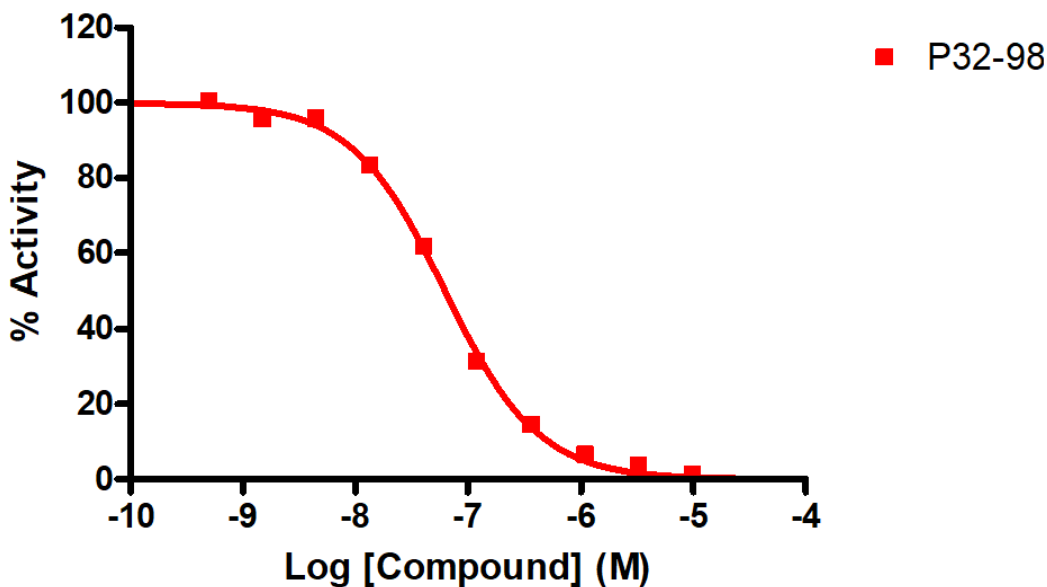
## Target Overview

CAT#	DPP-IV
Enzyme	Human Dipeptidyl peptidase IV
Synonym	DPP4, CD26, Adenosine deaminase complexing protein 2, ADCP2
Substrates	H-GP-AMC, 10 $\mu$ M

## Assay Properties

Readout	Enzymatic activity
Reference Compound IC50s	<b>Compounds IC50 (nM)</b> P32-98      120
Scientific Information	<a href="https://www.uniprot.org/uniprotkb/P27487/entry">P27487</a> (https://www.uniprot.org/uniprotkb/P27487/entry)
Screening Location	Malvern, PA, USA
Further information	More information can be found on our website <a href="#">Protease Assay Services</a> .

## Reference compound IC50 for DPP-IV



Let's Discover Together.

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

[https://www.reactionbiology.com/datasheet/dpp-iv\\_protea\\_malvern](https://www.reactionbiology.com/datasheet/dpp-iv_protea_malvern)

