

DPPIV drug discovery kit

The DPPIV Drug Discovery Kit is a complete assay system designed to screen DPPIV inhibitors, providing enough material to perform at least 96 assays. DPPIV (DPP4, CD26) is a member of the class of proteases known as prolyl peptidases, which cleave proteins after proline residues and is thought to play roles in diabetes, cancer, and autoimmune diseases, making it a target for drug discovery.

The kit contains both a chromogenic substrate (H-Gly-Pro-pNA; $K_m = 114 \mu\text{M}$) and a fluorogenic substrate (H-Gly-Pro-AMC; $K_m = 50 \mu\text{M}$).

Cleavage of the p-nitroaniline (pNA) from the colorimetric substrate increases absorbance at 405 nm. The fluorimetric assay is based on the cleavage of 7-amino-4-methylcoumarin (AMC) moiety from the C-terminus of the peptide substrate, which increases its fluorescence intensity at 460 nm. The kit is useful to screen inhibitors of DPPIV, a potential therapeutic target. A DPPIV inhibitor, P32/98 ($K_i = 130 \text{ nM}$), is included for use as a control. Other DPP enzymes are available for specificity profiling.

Citations: 24

[View Online »](#)

Ordering Information

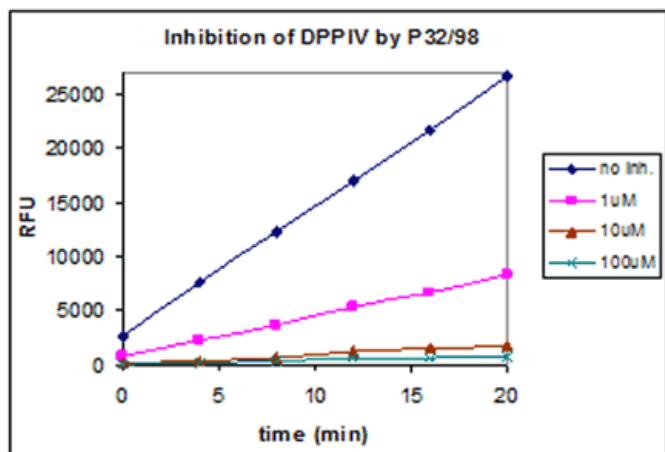
[Order Online »](#)

BML-AK499-0001	96 wells
----------------	----------

Manuals, SDS & CofA

[View Online »](#)

- Measures DPPIV activity in plasma, serum, urine, and saliva
- Allows correlation of DPPIV activity levels with disease states or drug treatments
- Allows determination of efficacy of DPPIV inhibitors



Plot data as Relative Fluorescence Units (RFU) versus time for each sample.

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	CD26, Dipeptidyl Peptidase IV, DPP4
Application	Activity assay, Colorimetric detection, Fluorescent detection, HTS
Compatibility	This product is compatible with the Absorbance 96 Plate Reader .
Contents	DPPIV enzyme (human, recombinant), pNA substrate, Calibration standard (p-nitroaniline), AMC substrate, Calibration standard (7-Amino-4-methylcoumarin), Inhibitor, Assay Buffer, 1/2-volume clear microplate
UniProt ID	P27487

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)